

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	A01	MmiRQP0923	MIMAT0000235	mmu-miR-202-3p	
QM043-01	A02	MmiRQP0421	MIMAT0000766	mmu-miR-335-5p	
QM043-01	A03	MmiRQP0029	MIMAT0000386	mmu-miR-106b-5p	
QM043-01	A04	MmiRQP0422	MIMAT0004704	mmu-miR-335-3p	
QM043-01	A05	MmiRQP0043	MIMAT0005858	mmu-miR-1197-3p	
QM043-01	A06	MmiRQP0426	MIMAT0000582	mmu-miR-338-3p	
QM043-01	A07	MmiRQP0044	MIMAT0000123	mmu-miR-1a-3p	
QM043-01	A08	MmiRQP0427	MIMAT0000584	mmu-miR-339-5p	
QM043-01	A09	MmiRQP0056	MIMAT0000246	mmu-miR-122-5p	
QM043-01	A10	MmiRQP0430	MIMAT0000667	mmu-miR-33-5p	
QM043-01	A11	MmiRQP0074	MIMAT0000134	mmu-miR-124-3p	
QM043-01	A12	MmiRQP0433	MIMAT0000586	mmu-miR-340-3p	
QM043-01	A13	MmiRQP0093	MIMAT0004528	mmu-miR-125a-3p	
QM043-01	A14	MmiRQP0434	MIMAT0004651	mmu-miR-340-5p	
QM043-01	A15	MmiRQP0096	MIMAT0000136	mmu-miR-125b-5p	
QM043-01	A16	MmiRQP0436	MIMAT0000590	mmu-miR-342-3p	
QM043-01	A17	MmiRQP0098	MIMAT0000137	mmu-miR-126-5p	
QM043-01	A18	MmiRQP0440	MIMAT0000542	mmu-miR-34a-5p	
QM043-01	A19	MmiRQP0099	MIMAT0000138	mmu-miR-126-3p	
QM043-01	A20	MmiRQP0446	MIMAT0000704	mmu-miR-361-5p	
QM043-01	A21	MmiRQP0111	MIMAT0000139	mmu-miR-127-3p	
QM043-01	A22	MmiRQP0449	MIMAT0000708	mmu-miR-363-3p	
QM043-01	A23	MmiRQP0125	MIMAT0000140	mmu-miR-128-3p	
QM043-01	A24	MmiRQP0451	MIMAT0000711	mmu-miR-365-3p	
QM043-01	B01	MmiRQP0895	MIMAT0000211	mmu-miR-182-5p	
QM043-01	B02	MmiRQP1074	MIMAT0005292	mmu-miR-582-3p	
QM043-01	B03	MmiRQP0897	MIMAT0007871	mmu-miR-1892	
QM043-01	B04	MmiRQP1075	MIMAT0005291	mmu-miR-582-5p	
QM043-01	B05	MmiRQP0898	MIMAT0007879	mmu-miR-1893	
QM043-01	B06	MmiRQP1076	MIMAT0003730	mmu-miR-592-5p	
QM043-01	B07	MmiRQP0899	MIMAT0007877	mmu-miR-1894-5p	
QM043-01	B08	MmiRQP1077	MIMAT0004942	mmu-miR-598-3p	
QM043-01	B09	MmiRQP0902	MIMAT0007873	mmu-miR-1896	
QM043-01	B10	MmiRQP1080	MIMAT0003737	mmu-miR-666-5p	
QM043-01	B11	MmiRQP0905	MIMAT0007875	mmu-miR-1898	
QM043-01	B12	MmiRQP1081	MIMAT0004823	mmu-miR-666-3p	
QM043-01	B13	MmiRQP0909	MIMAT0007870	mmu-miR-1900	
QM043-01	B14	MmiRQP1082	MIMAT0003734	mmu-miR-667-3p	
QM043-01	B15	MmiRQP0910	MIMAT0007880	mmu-miR-1901	
QM043-01	B16	MmiRQP1083	MIMAT0003732	mmu-miR-668-3p	
QM043-01	B17	MmiRQP0911	MIMAT0007863	mmu-miR-1902	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	B18	MmiRQP1088	MIMAT0005853	mmu-miR-669e-5p	
QM043-01	B19	MmiRQP0918	MIMAT0004859	mmu-miR-193b-3p	
QM043-01	B20	MmiRQP1089	MIMAT0005839	mmu-miR-669f-3p	
QM043-01	B21	MmiRQP0922	MIMAT0000234	mmu-miR-201-5p	
QM043-01	B22	MmiRQP1091	MIMAT0005841	mmu-miR-669h-5p	
QM043-01	B23	MmiRQP0924	MIMAT0004546	mmu-miR-202-5p	
QM043-01	B24	MmiRQP1096	MIMAT0003736	mmu-miR-670-5p	
QM043-01	C01	MmiRQP0137	MIMAT0000209	mmu-miR-129-5p	
QM043-01	C02	MmiRQP0454	MIMAT0003186	mmu-miR-369-3p	
QM043-01	C03	MmiRQP0139	MIMAT0000544	mmu-miR-129-2-3p	
QM043-01	C04	MmiRQP0455	MIMAT0003185	mmu-miR-369-5p	
QM043-01	C05	MmiRQP0161	MIMAT0000144	mmu-miR-132-3p	
QM043-01	C06	MmiRQP0456	MIMAT0001095	mmu-miR-370-3p	
QM043-01	C07	MmiRQP0168	MIMAT0000146	mmu-miR-134-5p	
QM043-01	C08	MmiRQP0464	MIMAT0003727	mmu-miR-374-5p	
QM043-01	C09	MmiRQP0175	MIMAT0000149	mmu-miR-137-3p	
QM043-01	C10	MmiRQP0466	MIMAT0000739	mmu-miR-375-3p	
QM043-01	C11	MmiRQP0180	MIMAT0000656	mmu-miR-139-5p	
QM043-01	C12	MmiRQP0472	MIMAT0000741	mmu-miR-377-3p	
QM043-01	C13	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM043-01	C14	MmiRQP0473	MIMAT0003151	mmu-miR-378-3p	
QM043-01	C15	MmiRQP0185	MIMAT0000154	mmu-miR-142-5p	
QM043-01	C16	MmiRQP0476	MIMAT0000743	mmu-miR-379-5p	
QM043-01	C17	MmiRQP0186	MIMAT0000155	mmu-miR-142-3p	
QM043-01	C18	MmiRQP0479	MIMAT0000746	mmu-miR-381-3p	
QM043-01	C19	MmiRQP0188	MIMAT0000247	mmu-miR-143-3p	
QM043-01	C20	MmiRQP0480	MIMAT0000747	mmu-miR-382-5p	
QM043-01	C21	MmiRQP0190	MIMAT0000156	mmu-miR-144-3p	
QM043-01	C22	MmiRQP0483	MIMAT0004746	mmu-miR-409-5p	
QM043-01	C23	MmiRQP0192	MIMAT0000157	mmu-miR-145-5p	
QM043-01	C24	MmiRQP0484	MIMAT0001090	mmu-miR-409-3p	
QM043-01	D01	MmiRQP0931	MIMAT0000659	mmu-miR-212-3p	
QM043-01	D02	MmiRQP1097	MIMAT0003735	mmu-miR-672-5p	
QM043-01	D03	MmiRQP0932	MIMAT0000904	mmu-miR-215-5p	
QM043-01	D04	MmiRQP1098	MIMAT0003739	mmu-miR-673-5p	
QM043-01	D05	MmiRQP0933	MIMAT0000679	mmu-miR-217-5p	
QM043-01	D06	MmiRQP1099	MIMAT0004824	mmu-miR-673-3p	
QM043-01	D07	MmiRQP0935	MIMAT0004863	mmu-miR-220	
QM043-01	D08	MmiRQP1104	MIMAT0003782	mmu-miR-676-3p	
QM043-01	D09	MmiRQP0936	MIMAT0000671	mmu-miR-224-5p	
QM043-01	D10	MmiRQP1105	MIMAT0003451	mmu-miR-677-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	D11	MmiRQP0939	MIMAT0004572	mmu-miR-290-3p	
QM043-01	D12	MmiRQP1111	MIMAT0003461	mmu-miR-683	
QM043-01	D13	MmiRQP0940	MIMAT0000366	mmu-miR-290-5p	
QM043-01	D14	MmiRQP1112	MIMAT0003462	mmu-miR-684	
QM043-01	D15	MmiRQP0942	MIMAT0003190	mmu-miR-291b-3p	
QM043-01	D16	MmiRQP1113	MIMAT0003463	mmu-miR-685	
QM043-01	D17	MmiRQP0944	MIMAT0000370	mmu-miR-292-3p	
QM043-01	D18	MmiRQP1114	MIMAT0003464	mmu-miR-686	
QM043-01	D19	MmiRQP0946	MIMAT0000371	mmu-miR-293-3p	
QM043-01	D20	MmiRQP1115	MIMAT0003466	mmu-miR-687	
QM043-01	D21	MmiRQP0947	MIMAT0000372	mmu-miR-294-3p	
QM043-01	D22	MmiRQP1116	MIMAT0003467	mmu-miR-688	
QM043-01	D23	MmiRQP0949	MIMAT0000373	mmu-miR-295-3p	
QM043-01	D24	MmiRQP1117	MIMAT0003468	mmu-miR-689	
QM043-01	E01	MmiRQP0202	MIMAT0004857	mmu-miR-147-3p	
QM043-01	E02	MmiRQP0485	MIMAT0001091	mmu-miR-410-3p	
QM043-01	E03	MmiRQP0208	MIMAT0000159	mmu-miR-149-5p	
QM043-01	E04	MmiRQP0486	MIMAT0004747	mmu-miR-411-5p	
QM043-01	E05	MmiRQP0210	MIMAT0000160	mmu-miR-150-5p	
QM043-01	E06	MmiRQP0489	MIMAT0004869	mmu-miR-421-3p	
QM043-01	E07	MmiRQP0212	MIMAT0004536	mmu-miR-151-5p	
QM043-01	E08	MmiRQP0491	MIMAT0003454	mmu-miR-423-3p	
QM043-01	E09	MmiRQP0214	MIMAT0000163	mmu-miR-153-3p	
QM043-01	E10	MmiRQP0492	MIMAT0004825	mmu-miR-423-5p	
QM043-01	E11	MmiRQP0219	MIMAT0000164	mmu-miR-154-5p	
QM043-01	E12	MmiRQP0495	MIMAT0004750	mmu-miR-425-5p	
QM043-01	E13	MmiRQP0223	MIMAT0000526	mmu-miR-15a-5p	
QM043-01	E14	MmiRQP0499	MIMAT0001418	mmu-miR-431-5p	
QM043-01	E15	MmiRQP0225	MIMAT0000124	mmu-miR-15b-5p	
QM043-01	E16	MmiRQP0502	MIMAT0001420	mmu-miR-433-3p	
QM043-01	E17	MmiRQP0227	MIMAT0000527	mmu-miR-16-5p	
QM043-01	E18	MmiRQP0503	MIMAT0001533	mmu-miR-448-3p	
QM043-01	E19	MmiRQP0244	MIMAT0000212	mmu-miR-183-5p	
QM043-01	E20	MmiRQP0504	MIMAT0001542	mmu-miR-449a-5p	
QM043-01	E21	MmiRQP0245	MIMAT0000213	mmu-miR-184-3p	
QM043-01	E22	MmiRQP0506	MIMAT0001546	mmu-miR-450a-5p	
QM043-01	E23	MmiRQP0247	MIMAT0000214	mmu-miR-185-5p	
QM043-01	E24	MmiRQP0509	MIMAT0001632	mmu-miR-451	
QM043-01	F01	MmiRQP0954	MIMAT0004827	mmu-miR-297b-3p	mmu-miR-297c-3p mmu-miR-297a-3p
QM043-01	F02	MmiRQP1119	MIMAT0003470	mmu-miR-691	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	F03	MmiRQP0955	MIMAT0000376	mmu-miR-298-5p	
QM043-01	F04	MmiRQP1122	MIMAT0004189	mmu-miR-693-3p	
QM043-01	F05	MmiRQP0956	MIMAT0004577	mmu-miR-299-3p	
QM043-01	F06	MmiRQP1123	MIMAT0003474	mmu-miR-694	
QM043-01	F07	MmiRQP0958	MIMAT0000378	mmu-miR-300-3p	
QM043-01	F08	MmiRQP1124	MIMAT0003481	mmu-miR-695	
QM043-01	F09	MmiRQP0967	MIMAT0000556	mmu-miR-324-3p	
QM043-01	F10	MmiRQP1125	MIMAT0003483	mmu-miR-696	
QM043-01	F11	MmiRQP0968	MIMAT0004640	mmu-miR-325-3p	
QM043-01	F12	MmiRQP1126	MIMAT0003487	mmu-miR-697	
QM043-01	F13	MmiRQP0969	MIMAT0004867	mmu-miR-327	
QM043-01	F14	MmiRQP1127	MIMAT0003488	mmu-miR-698	
QM043-01	F15	MmiRQP0971	MIMAT0004644	mmu-miR-337-5p	
QM043-01	F16	MmiRQP1129	MIMAT0003490	mmu-miR-700-3p	
QM043-01	F17	MmiRQP0972	MIMAT0000578	mmu-miR-337-3p	
QM043-01	F18	MmiRQP1130	MIMAT0003491	mmu-miR-701-5p	
QM043-01	F19	MmiRQP0973	MIMAT0004649	mmu-miR-339-3p	
QM043-01	F20	MmiRQP1132	MIMAT0003493	mmu-miR-703	
QM043-01	F21	MmiRQP0974	MIMAT0004868	mmu-miR-343	
QM043-01	F22	MmiRQP1134	MIMAT0003495	mmu-miR-705	
QM043-01	F23	MmiRQP0975	MIMAT0000593	mmu-miR-344-3p	
QM043-01	F24	MmiRQP1135	MIMAT0003496	mmu-miR-706	
QM043-01	G01	MmiRQP0248	MIMAT0000215	mmu-miR-186-5p	
QM043-01	G02	MmiRQP0519	MIMAT0003127	mmu-miR-484	
QM043-01	G03	MmiRQP0252	MIMAT0000217	mmu-miR-188-5p	
QM043-01	G04	MmiRQP0520	MIMAT0003128	mmu-miR-485-5p	
QM043-01	G05	MmiRQP0253	MIMAT0004541	mmu-miR-188-3p	
QM043-01	G06	MmiRQP0525	MIMAT0003184	mmu-miR-487b-3p	
QM043-01	G07	MmiRQP0258	MIMAT0000220	mmu-miR-190-5p	
QM043-01	G08	MmiRQP0529	MIMAT0003780	mmu-miR-490-3p	
QM043-01	G09	MmiRQP0262	MIMAT0004852	mmu-miR-190b-5p	
QM043-01	G10	MmiRQP0531	MIMAT0003486	mmu-miR-491-5p	
QM043-01	G11	MmiRQP0263	MIMAT0000221	mmu-miR-191-5p	
QM043-01	G12	MmiRQP0536	MIMAT0003182	mmu-miR-494-3p	
QM043-01	G13	MmiRQP0274	MIMAT0000517	mmu-miR-192-5p	
QM043-01	G14	MmiRQP0537	MIMAT0003456	mmu-miR-495-3p	
QM043-01	G15	MmiRQP0276	MIMAT0000223	mmu-miR-193-3p	
QM043-01	G16	MmiRQP0538	MIMAT0003738	mmu-miR-496-3p	
QM043-01	G17	MmiRQP0280	MIMAT0000224	mmu-miR-194-5p	
QM043-01	G18	MmiRQP0543	MIMAT0003482	mmu-miR-499-5p	
QM043-01	G19	MmiRQP0283	MIMAT0000225	mmu-miR-195-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	G20	MmiRQP0551	MIMAT0004889	mmu-miR-504-5p	
QM043-01	G21	MmiRQP0290	MIMAT0000229	mmu-miR-199a-5p	
QM043-01	G22	MmiRQP0615	MIMAT0002889	mmu-miR-532-5p	
QM043-01	G23	MmiRQP0300	MIMAT0000233	mmu-miR-200b-3p	
QM043-01	G24	MmiRQP0616	MIMAT0004781	mmu-miR-532-3p	
QM043-01	H01	MmiRQP0977	MIMAT0000595	mmu-miR-345-5p	
QM043-01	H02	MmiRQP1136	MIMAT0003497	mmu-miR-707	
QM043-01	H03	MmiRQP0978	MIMAT0000597	mmu-miR-346-5p	
QM043-01	H04	MmiRQP1140	MIMAT0003502	mmu-miR-712-5p	
QM043-01	H05	MmiRQP0979	MIMAT0004581	mmu-miR-34b-3p	
QM043-01	H06	MmiRQP1143	MIMAT0003505	mmu-miR-714	
QM043-01	H07	MmiRQP0982	MIMAT0000605	mmu-miR-350-3p	
QM043-01	H08	MmiRQP1144	MIMAT0003506	mmu-miR-715	
QM043-01	H09	MmiRQP0984	MIMAT0004684	mmu-miR-362-3p	
QM043-01	H10	MmiRQP1145	MIMAT0003510	mmu-miR-717	
QM043-01	H11	MmiRQP0985	MIMAT0000706	mmu-miR-362-5p	
QM043-01	H12	MmiRQP1146	MIMAT0003514	mmu-miR-718	
QM043-01	H13	MmiRQP0987	MIMAT0000740	mmu-miR-376a-3p	
QM043-01	H14	MmiRQP1147	MIMAT0003465	mmu-miR-719	
QM043-01	H15	MmiRQP0989	MIMAT0001092	mmu-miR-376b-3p	
QM043-01	H16	MmiRQP1148	MIMAT0003484	mmu-miR-720	
QM043-01	H17	MmiRQP0990	MIMAT0003183	mmu-miR-376c-3p	
QM043-01	H18	MmiRQP1149	MIMAT0003515	mmu-miR-721	
QM043-01	H19	MmiRQP0992	MIMAT0000744	mmu-miR-380-5p	
QM043-01	H20	MmiRQP1150	MIMAT0004236	mmu-miR-741-3p	
QM043-01	H21	MmiRQP0993	MIMAT0000745	mmu-miR-380-3p	
QM043-01	H22	MmiRQP1151	MIMAT0004237	mmu-miR-742-3p	
QM043-01	H23	MmiRQP0995	MIMAT0000748	mmu-miR-383-5p	
QM043-01	H24	MmiRQP1153	MIMAT0004238	mmu-miR-743a-3p	
QM043-01	I01	MmiRQP0302	MIMAT0000657	mmu-miR-200c-3p	
QM043-01	I02	MmiRQP0617	MIMAT0003169	mmu-miR-539-5p	
QM043-01	I03	MmiRQP0305	MIMAT0000236	mmu-miR-203-3p	
QM043-01	I04	MmiRQP0621	MIMAT0003172	mmu-miR-542-3p	
QM043-01	I05	MmiRQP0307	MIMAT0000238	mmu-miR-205-5p	
QM043-01	I06	MmiRQP0622	MIMAT0003168	mmu-miR-543-3p	
QM043-01	I07	MmiRQP0308	MIMAT0000239	mmu-miR-206-3p	
QM043-01	I08	MmiRQP0651	MIMAT0003890	mmu-miR-551b-3p	
QM043-01	I09	MmiRQP0309	MIMAT0000520	mmu-miR-208a-3p	
QM043-01	I10	MmiRQP0667	MIMAT0004892	mmu-miR-568	
QM043-01	I11	MmiRQP0310	MIMAT0004939	mmu-miR-208b-3p	
QM043-01	I12	MmiRQP0673	MIMAT0004894	mmu-miR-574-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	I13	MmiRQP0316	MIMAT0000530	mmu-miR-21-5p	
QM043-01	I14	MmiRQP0674	MIMAT0004893	mmu-miR-574-5p	
QM043-01	I15	MmiRQP0317	MIMAT0000658	mmu-miR-210-3p	
QM043-01	I16	MmiRQP0693	MIMAT0004895	mmu-miR-590-5p	
QM043-01	I17	MmiRQP0320	MIMAT0000661	mmu-miR-214-3p	
QM043-01	I18	MmiRQP0694	MIMAT0004896	mmu-miR-590-3p	
QM043-01	I19	MmiRQP0324	MIMAT0003729	mmu-miR-216b-5p	
QM043-01	I20	MmiRQP0719	MIMAT0004837	mmu-miR-615-5p	
QM043-01	I21	MmiRQP0327	MIMAT0000663	mmu-miR-218-5p	
QM043-01	I22	MmiRQP0720	MIMAT0003783	mmu-miR-615-3p	
QM043-01	I23	MmiRQP0330	MIMAT0000664	mmu-miR-219-5p	
QM043-01	I24	MmiRQP0762	MIMAT0003711	mmu-miR-652-3p	
QM043-01	J01	MmiRQP0996	MIMAT0001076	mmu-miR-384-3p	
QM043-01	J02	MmiRQP1154	MIMAT0004840	mmu-miR-743b-3p	
QM043-01	J03	MmiRQP0997	MIMAT0001094	mmu-miR-412-3p	
QM043-01	J04	MmiRQP1155	MIMAT0004839	mmu-miR-743b-5p	
QM043-01	J05	MmiRQP0999	MIMAT0001537	mmu-miR-429-3p	
QM043-01	J06	MmiRQP1156	MIMAT0003889	mmu-miR-758-3p	
QM043-01	J07	MmiRQP1001	MIMAT0001421	mmu-miR-434-5p	
QM043-01	J08	MmiRQP1157	MIMAT0003897	mmu-miR-759	
QM043-01	J09	MmiRQP1002	MIMAT0001422	mmu-miR-434-3p	
QM043-01	J10	MmiRQP1158	MIMAT0003893	mmu-miR-761	
QM043-01	J11	MmiRQP1006	MIMAT0003512	mmu-miR-450b-3p	
QM043-01	J12	MmiRQP1160	MIMAT0003896	mmu-miR-763	
QM043-01	J13	MmiRQP1007	MIMAT0003511	mmu-miR-450b-5p	
QM043-01	J14	MmiRQP1161	MIMAT0003895	mmu-miR-764-3p	
QM043-01	J15	MmiRQP1010	MIMAT0003742	mmu-miR-455-3p	
QM043-01	J16	MmiRQP1162	MIMAT0003894	mmu-miR-764-5p	
QM043-01	J17	MmiRQP1012	MIMAT0004758	mmu-miR-463-3p	
QM043-01	J18	MmiRQP1163	MIMAT0004822	mmu-miR-770-5p	
QM043-01	J19	MmiRQP1013	MIMAT0002105	mmu-miR-464	
QM043-01	J20	MmiRQP1165	MIMAT0004188	mmu-miR-802-5p	
QM043-01	J21	MmiRQP1014	MIMAT0002106	mmu-miR-465a-5p	
QM043-01	J22	MmiRQP1166	MIMAT0004210	mmu-miR-804	
QM043-01	J23	MmiRQP1023	MIMAT0004930	mmu-miR-466d-5p	
QM043-01	J24	MmiRQP1167	MIMAT0004211	mmu-miR-805	
QM043-01	K01	MmiRQP0332	MIMAT0000531	mmu-miR-22-3p	
QM043-01	K02	MmiRQP0763	MIMAT0004943	mmu-miR-653-5p	
QM043-01	K03	MmiRQP0338	MIMAT0000669	mmu-miR-221-3p	
QM043-01	K04	MmiRQP0764	MIMAT0004898	mmu-miR-654-3p	
QM043-01	K05	MmiRQP0339	MIMAT0000670	mmu-miR-222-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	K06	MmiRQP0781	MIMAT0004821	mmu-miR-671-3p	
QM043-01	K07	MmiRQP0342	MIMAT0000665	mmu-miR-223-3p	
QM043-01	K08	MmiRQP0783	MIMAT0004828	mmu-miR-708-5p	
QM043-01	K09	MmiRQP0349	MIMAT0000219	mmu-miR-24-3p	
QM043-01	K10	MmiRQP0790	MIMAT0004187	mmu-miR-744-5p	
QM043-01	K11	MmiRQP0352	MIMAT0000652	mmu-miR-25-3p	
QM043-01	K12	MmiRQP0792	MIMAT0003898	mmu-miR-760-3p	
QM043-01	K13	MmiRQP0354	MIMAT0000533	mmu-miR-26a-5p	
QM043-01	K14	MmiRQP0803	MIMAT0004936	mmu-miR-873-5p	
QM043-01	K15	MmiRQP0357	MIMAT0000534	mmu-miR-26b-5p	
QM043-01	K16	MmiRQP0804	MIMAT0004853	mmu-miR-874-3p	
QM043-01	K17	MmiRQP0362	MIMAT0000653	mmu-miR-28-5p	
QM043-01	K18	MmiRQP0806	MIMAT0004937	mmu-miR-875-5p	
QM043-01	K19	MmiRQP0364	MIMAT0000374	mmu-miR-296-5p	
QM043-01	K20	MmiRQP0809	MIMAT0004861	mmu-miR-877-5p	
QM043-01	K21	MmiRQP0365	MIMAT0004576	mmu-miR-296-3p	
QM043-01	K22	MmiRQP0825	MIMAT0000142	mmu-miR-9-5p	
QM043-01	K23	MmiRQP0373	MIMAT0000127	mmu-miR-29b-3p	
QM043-01	K24	MmiRQP0837	MIMAT0000540	mmu-miR-93-5p	
QM043-01	L01	MmiRQP1026	MIMAT0004882	mmu-miR-466f-3p	
QM043-01	L02	MmiRQP1168	MIMAT0004841	mmu-miR-871-5p	
QM043-01	L03	MmiRQP1029	MIMAT0004883	mmu-miR-466g	
QM043-01	L04	MmiRQP1170	MIMAT0004938	mmu-miR-875-3p	
QM043-01	L05	MmiRQP1031	MIMAT0005834	mmu-miR-466i-3p	
QM043-01	L06	MmiRQP1171	MIMAT0004855	mmu-miR-876-3p	
QM043-01	L07	MmiRQP1033	MIMAT0005845	mmu-miR-466k	
QM043-01	L08	MmiRQP1172	MIMAT0004854	mmu-miR-876-5p	
QM043-01	L09	MmiRQP1034	MIMAT0005830	mmu-miR-466l-3p	
QM043-01	L10	MmiRQP1174	MIMAT0004933	mmu-miR-878-3p	
QM043-01	L11	MmiRQP1037	MIMAT0005448	mmu-miR-467b-5p	
QM043-01	L12	MmiRQP1175	MIMAT0004932	mmu-miR-878-5p	
QM043-01	L13	MmiRQP1043	MIMAT0005846	mmu-miR-467f	
QM043-01	L14	MmiRQP1179	MIMAT0004846	mmu-miR-881-3p	
QM043-01	L15	MmiRQP1047	MIMAT0002110	mmu-miR-469	
QM043-01	L16	MmiRQP1181	MIMAT0004847	mmu-miR-882	
QM043-01	L17	MmiRQP1049	MIMAT0002111	mmu-miR-470-5p	
QM043-01	L18	MmiRQP1183	MIMAT0004848	mmu-miR-883a-5p	
QM043-01	L19	MmiRQP1050	MIMAT0002112	mmu-miR-471-5p	
QM043-01	L20	MmiRQP1224	MIMAT0000538	mmu-miR-31-5p	
QM043-01	L21	MmiRQP1055	MIMAT0003450	mmu-miR-488-3p	
QM043-01	L22	MmiRQP1226	MIMAT0000548	mmu-miR-322-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	L23	MmiRQP1056	MIMAT0003112	mmu-miR-489-3p	
QM043-01	L24	MmiRQP1229	MIMAT0000559	mmu-miR-326-3p	
QM043-01	M01	MmiRQP0381	MIMAT0000380	mmu-miR-302a-3p	
QM043-01	M02	MmiRQP0852	MIMAT0000541	mmu-miR-96-5p	
QM043-01	M03	MmiRQP0393	MIMAT0000130	mmu-miR-30b-5p	
QM043-01	M04	MmiRQP0857	MIMAT0000132	mmu-miR-99b-5p	
QM043-01	M05	MmiRQP0396	MIMAT0000514	mmu-miR-30c-5p	
QM043-01	M06	MmiRQP0861	MIMAT0004856	mmu-miR-105	
QM043-01	M07	MmiRQP0404	MIMAT0000654	mmu-miR-32-5p	
QM043-01	M08	MmiRQP0868	MIMAT0005849	mmu-miR-1191	
QM043-01	M09	MmiRQP0405	MIMAT0000666	mmu-miR-320-3p	
QM043-01	M10	MmiRQP0869	MIMAT0005850	mmu-miR-1192	
QM043-01	M11	MmiRQP0409	MIMAT0004638	mmu-miR-323-5p	
QM043-01	M12	MmiRQP0870	MIMAT0005851	mmu-miR-1193-3p	
QM043-01	M13	MmiRQP0410	MIMAT0000551	mmu-miR-323-3p	
QM043-01	M14	MmiRQP0871	MIMAT0005852	mmu-miR-1194	
QM043-01	M15	MmiRQP0412	MIMAT0000555	mmu-miR-324-5p	
QM043-01	M16	MmiRQP0875	MIMAT0005860	mmu-miR-1199-5p	
QM043-01	M17	MmiRQP0415	MIMAT0000565	mmu-miR-328-3p	
QM043-01	M18	MmiRQP0878	MIMAT0005835	mmu-miR-1b-5p	
QM043-01	M19	MmiRQP0418	MIMAT0004642	mmu-miR-330-5p	
QM043-01	M20	MmiRQP0881	MIMAT0000148	mmu-miR-136-5p	
QM043-01	M21	MmiRQP0419	MIMAT0004643	mmu-miR-331-5p	
QM043-01	M22	MmiRQP0889	MIMAT0000161	mmu-miR-151-3p	
QM043-01	M23	MmiRQP0420	MIMAT0000571	mmu-miR-331-3p	
QM043-01	M24	MmiRQP0890	MIMAT0000165	mmu-miR-155-5p	
QM043-01	N01	MmiRQP1058	MIMAT0003453	mmu-miR-497-5p	
QM043-01	N02	MmiRQP1231	MIMAT0000567	mmu-miR-329-3p	
QM043-01	N03	MmiRQP1059	MIMAT0003507	mmu-miR-500-3p	
QM043-01	N04	MmiRQP1238	MIMAT0000588	mmu-miR-341-3p	
QM043-01	N05	MmiRQP1060	MIMAT0003508	mmu-miR-501-5p	
QM043-01	N06	MmiRQP1239	MIMAT0004653	mmu-miR-342-5p	
QM043-01	N07	MmiRQP1061	MIMAT0003509	mmu-miR-501-3p	
QM043-01	N08	MmiRQP1259	MIMAT0004745	mmu-miR-384-5p	
QM043-01	N09	MmiRQP1063	MIMAT0003188	mmu-miR-503-5p	
QM043-01	N10	MmiRQP1279	MIMAT0012774	mmu-miR-664-3p	
QM043-01	N11	MmiRQP1064	MIMAT0003513	mmu-miR-505-3p	
QM043-01	N12	MmiRQP1281	MIMAT0003740	mmu-miR-674-5p	
QM043-01	N13	MmiRQP1065	MIMAT0004890	mmu-miR-509-5p	
QM043-01	N14	MmiRQP1289	MIMAT0004934	mmu-miR-872-5p	
QM043-01	N15	MmiRQP1068	MIMAT0003167	mmu-miR-540-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	N16	MmiRQP1335	MIMAT0009391	mmu-miR-1928	
QM043-01	N17	MmiRQP1069	MIMAT0004786	mmu-miR-540-5p	
QM043-01	N18	MmiRQP1337	MIMAT0009393	mmu-miR-1930-5p	
QM043-01	N19	MmiRQP1070	MIMAT0003170	mmu-miR-541-5p	
QM043-01	N20	MmiRQP1339	MIMAT0009395	mmu-miR-1932	
QM043-01	N21	MmiRQP1071	MIMAT0004941	mmu-miR-544-3p	
QM043-01	N22	MmiRQP1340	MIMAT0009396	mmu-miR-1933-5p	
QM043-01	N23	MmiRQP1073	MIMAT0003173	mmu-miR-547-3p	
QM043-01	N24	MmiRQP1346	MIMAT0009403	mmu-miR-1939	
QM043-01	O01	MmiRQP2556	MIMAT0020636	mmu-miR-3572	
QM043-01	O02	MmiRQP2673	MIMAT0022836	mmu-miR-299b-5p	
QM043-01	O03	MmiRQP2557	MIMAT0020637	mmu-miR-5126	
QM043-01	O04	MmiRQP2674	MIMAT0022837	mmu-miR-299b-3p	
QM043-01	O05	MmiRQP2558	MIMAT0020638	mmu-miR-5127	
QM043-01	O06	MmiRQP2675	MIMAT0022355	mmu-miR-5615-5p	
QM043-01	O07	MmiRQP2559	MIMAT0020639	mmu-miR-5128	
QM043-01	O08	MmiRQP2676	MIMAT0022356	mmu-miR-5615-3p	
QM043-01	O09	MmiRQP2560	MIMAT0020640	mmu-miR-5129	
QM043-01	O10	MmiRQP2677	MIMAT0022357	mmu-miR-1231-5p	
QM043-01	O11	MmiRQP2561	MIMAT0020641	mmu-miR-5130	
QM043-01	O12	MmiRQP2678	MIMAT0022358	mmu-miR-1231-3p	
QM043-01	O13	MmiRQP2562	MIMAT0020642	mmu-miR-5131	
QM043-01	O14	MmiRQP2679	MIMAT0022359	mmu-miR-5616-5p	
QM043-01	O15	MmiRQP2563	MIMAT0020643	mmu-miR-5132	
QM043-01	O16	MmiRQP2680	MIMAT0022360	mmu-miR-5616-3p	
QM043-01	O17	MmiRQP2564	MIMAT0020644	mmu-miR-5133	
QM043-01	O18	MmiRQP2681	MIMAT0022361	mmu-miR-5617-5p	
QM043-01	O19	MmiRQP2566	MIMAT0020646	mmu-miR-5135	
QM043-01	O20	MmiRQP2682	MIMAT0022362	mmu-miR-5617-3p	
QM043-01	O21	MmiRQP2671	MIMAT0022353	mmu-miR-3544-5p	
QM043-01	O22	MmiRQP2683	MIMAT0022363	mmu-miR-5618-5p	
QM043-01	O23	MmiRQP2672	MIMAT0022354	mmu-miR-3544-3p	
QM043-01	O24	MmiRQP2684	MIMAT0022364	mmu-miR-5618-3p	
QM043-01	P01	NC	NC	NC	
QM043-01	P02	NC	NC	NC	
QM043-01	P03	NC	NC	NC	
QM043-01	P04	NC	NC	NC	
QM043-01	P05	MmiRQP9002	NR_003027	MK1	
QM043-01	P06	MmiRQP9002	NR_003027	MK1	
QM043-01	P07	MmiRQP9012	NR_028128	MK2	
QM043-01	P08	MmiRQP9012	NR_028128	MK2	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-01	P09	MmiRQP9022	NR_028554	MK3	
QM043-01	P10	MmiRQP9022	NR_028554	MK3	
QM043-01	P11	MmiRQP9042	NR_028550	MK4	
QM043-01	P12	MmiRQP9042	NR_028550	MK4	
QM043-01	P13	MmiRQP9052	NR_028541	MK5	
QM043-01	P14	MmiRQP9052	NR_028541	MK5	
QM043-01	P15	MmiRQP9062	NR_028091	MK6	
QM043-01	P16	MmiRQP9062	NR_028091	MK6	
QM043-01	P17	RT	RT	RT	
QM043-01	P18	RT	RT	RT	
QM043-01	P19	RT	RT	RT	
QM043-01	P20	RT	RT	RT	
QM043-01	P21	PCR	PCR	PCR	
QM043-01	P22	PCR	PCR	PCR	
QM043-01	P23	PCR	PCR	PCR	
QM043-01	P24	PCR	PCR	PCR	
QM043-02	A01	MmiRQP1347	MIMAT0009404	mmu-miR-1940	
QM043-02	A02	MmiRQP0166	MIMAT0000145	mmu-miR-133a-3p	
QM043-02	A03	MmiRQP1348	MIMAT0009405	mmu-miR-1941-5p	
QM043-02	A04	MmiRQP0167	MIMAT0000769	mmu-miR-133b-3p	
QM043-02	A05	MmiRQP1349	MIMAT0009407	mmu-miR-1942	
QM043-02	A06	MmiRQP0170	MIMAT0000147	mmu-miR-135a-5p	
QM043-02	A07	MmiRQP1351	MIMAT0009409	mmu-miR-1944	
QM043-02	A08	MmiRQP0172	MIMAT0000612	mmu-miR-135b-5p	
QM043-02	A09	MmiRQP1352	MIMAT0009410	mmu-miR-1945	
QM043-02	A10	MmiRQP0230	MIMAT0000649	mmu-miR-17-5p	
QM043-02	A11	MmiRQP1357	MIMAT0009416	mmu-miR-1949	
QM043-02	A12	MmiRQP0232	MIMAT0000210	mmu-miR-181a-5p	
QM043-02	A13	MmiRQP1358	MIMAT0009417	mmu-miR-1950	
QM043-02	A14	MmiRQP0234	MIMAT0000673	mmu-miR-181b-5p	
QM043-02	A15	MmiRQP1360	MIMAT0009419	mmu-miR-669m-3p	
QM043-02	A16	MmiRQP0235	MIMAT0000674	mmu-miR-181c-5p	
QM043-02	A17	MmiRQP1362	MIMAT0009422	mmu-miR-1951	
QM043-02	A18	MmiRQP0237	MIMAT0004324	mmu-miR-181d-5p	
QM043-02	A19	MmiRQP1363	MIMAT0009423	mmu-miR-1952	
QM043-02	A20	MmiRQP0284	MIMAT0000518	mmu-miR-196a-5p	
QM043-02	A21	MmiRQP1365	MIMAT0009425	mmu-miR-1954	
QM043-02	A22	MmiRQP0293	MIMAT0000651	mmu-miR-19a-3p	
QM043-02	A23	MmiRQP1366	MIMAT0009426	mmu-miR-1955-5p	
QM043-02	A24	MmiRQP0295	MIMAT0000513	mmu-miR-19b-3p	
QM043-02	B01	MmiRQP2744	MIMAT0017064	mmu-miR-135a-2-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	B02	MmiRQP1039	MIMAT0004887	mmu-miR-467d-3p	mmu-miR-467a-3p
QM043-02	B03	MmiRQP1725	MIMAT0004877	mmu-miR-466c-5p	
QM043-02	B04	MmiRQP1042	MIMAT0005294	mmu-miR-467e-3p	
QM043-02	B05	MmiRQP1723	MIMAT0003477	mmu-miR-669a-5p	mmu-miR-669p-5p
QM043-02	B06	MmiRQP1048	MIMAT0004760	mmu-miR-470-3p	
QM043-02	B07	MmiRQP1415	MIMAT0009461	mmu-miR-1982.2	
QM043-02	B08	MmiRQP1052	MIMAT0003120	mmu-miR-483-3p	
QM043-02	B09	MmiRQP1726	MIMAT0004878	mmu-miR-466c-3p	mmu-miR-466e-3p mmu-miR-466b-3p mmu-miR-466a-3p
QM043-02	B10	MmiRQP1054	MIMAT0003449	mmu-miR-488-5p	
QM043-02	B11	MmiRQP2569	MIMAT0017245	mmu-miR-760-5p	
QM043-02	B12	MmiRQP1100	MIMAT0003741	mmu-miR-674-3p	
QM043-02	B13	MmiRQP0003	MIMAT0004621	mmu-let-7b-3p	
QM043-02	B14	MmiRQP1103	MIMAT0003781	mmu-miR-676-5p	
QM043-02	B15	MmiRQP0008	MIMAT0000384	mmu-let-7d-3p	
QM043-02	B16	MmiRQP1128	MIMAT0003489	mmu-miR-699	
QM043-02	B17	MmiRQP0011	MIMAT0004623	mmu-let-7f-1-3p	
QM043-02	B18	MmiRQP1141	MIMAT0003743	mmu-miR-712-3p	
QM043-02	B19	MmiRQP0016	MIMAT0004520	mmu-let-7i-3p	
QM043-02	B20	MmiRQP1152	MIMAT0004838	mmu-miR-742-5p	
QM043-02	B21	MmiRQP0028	MIMAT0004582	mmu-miR-106b-3p	
QM043-02	B22	MmiRQP1169	MIMAT0004935	mmu-miR-872-3p	
QM043-02	B23	MmiRQP0031	MIMAT0004659	mmu-miR-10a-3p	
QM043-02	B24	MmiRQP1173	MIMAT0004862	mmu-miR-877-3p	
QM043-02	C01	MmiRQP1368	MIMAT0009428	mmu-miR-1956	
QM043-02	C02	MmiRQP0306	MIMAT0000237	mmu-miR-204-5p	
QM043-02	C03	MmiRQP1370	MIMAT0009430	mmu-miR-1957	
QM043-02	C04	MmiRQP0344	MIMAT0000532	mmu-miR-23a-3p	
QM043-02	C05	MmiRQP1371	MIMAT0009431	mmu-miR-1958	
QM043-02	C06	MmiRQP0346	MIMAT0000125	mmu-miR-23b-3p	
QM043-02	C07	MmiRQP1372	MIMAT0009432	mmu-miR-1959	
QM043-02	C08	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM043-02	C09	MmiRQP1374	MIMAT0009434	mmu-miR-1961	
QM043-02	C10	MmiRQP0361	MIMAT0000126	mmu-miR-27b-3p	
QM043-02	C11	MmiRQP1376	MIMAT0009436	mmu-miR-1963	
QM043-02	C12	MmiRQP0371	MIMAT0000535	mmu-miR-29a-3p	
QM043-02	C13	MmiRQP1378	MIMAT0009438	mmu-miR-1965	
QM043-02	C14	MmiRQP0375	MIMAT0000536	mmu-miR-29c-3p	
QM043-02	C15	MmiRQP1379	MIMAT0009439	mmu-miR-1966	
QM043-02	C16	MmiRQP0378	MIMAT0000379	mmu-miR-301a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	C17	MmiRQP1381	MIMAT0009441	mmu-miR-1968-5p	
QM043-02	C18	MmiRQP0391	MIMAT0000128	mmu-miR-30a-5p	
QM043-02	C19	MmiRQP1382	MIMAT0009442	mmu-miR-1969	
QM043-02	C20	MmiRQP0398	MIMAT0000515	mmu-miR-30d-5p	
QM043-02	C21	MmiRQP1384	MIMAT0009444	mmu-miR-1970	
QM043-02	C22	MmiRQP0400	MIMAT0000248	mmu-miR-30e-5p	
QM043-02	C23	MmiRQP1385	MIMAT0009445	mmu-miR-1274a	
QM043-02	C24	MmiRQP0786	MIMAT0000677	mmu-miR-7a-5p	
QM043-02	D01	MmiRQP0073	MIMAT0004527	mmu-miR-124-5p	
QM043-02	D02	MmiRQP1177	MIMAT0004843	mmu-miR-879-3p	
QM043-02	D03	MmiRQP0079	MIMAT0014800	mmu-miR-1247-5p	
QM043-02	D04	MmiRQP1199	MIMAT0004544	mmu-miR-193-5p	
QM043-02	D05	MmiRQP0081	MIMAT0010560	mmu-miR-1249-3p	
QM043-02	D06	MmiRQP1214	MIMAT0005440	mmu-miR-24-2-5p	
QM043-02	D07	MmiRQP0145	MIMAT0014809	mmu-miR-1298-5p	
QM043-02	D08	MmiRQP1228	MIMAT0000558	mmu-miR-325-5p	
QM043-02	D09	MmiRQP0169	MIMAT0004531	mmu-miR-135a-1-3p	
QM043-02	D10	MmiRQP1390	MIMAT0009459	mmu-miR-1982-5p	
QM043-02	D11	MmiRQP0182	MIMAT0000152	mmu-miR-140-3p	
QM043-02	D12	MmiRQP1623	MIMAT0015219	mmu-miR-3475	
QM043-02	D13	MmiRQP0203	MIMAT0004617	mmu-miR-148a-5p	
QM043-02	D14	MmiRQP1624	MIMAT0014815	mmu-miR-3099-5p	
QM043-02	D15	MmiRQP0224	MIMAT0004521	mmu-miR-15b-3p	
QM043-02	D16	MmiRQP1625	MIMAT0013803	mmu-miR-2861	
QM043-02	D17	MmiRQP0231	MIMAT0000660	mmu-miR-181a-1-3p	
QM043-02	D18	MmiRQP1626	MIMAT0014852	mmu-miR-3072-5p	
QM043-02	D19	MmiRQP0243	MIMAT0004539	mmu-miR-183-3p	
QM043-02	D20	MmiRQP1628	MIMAT0015641	mmu-miR-3470b	
QM043-02	D21	MmiRQP0297	MIMAT0004619	mmu-miR-200a-5p	
QM043-02	D22	MmiRQP1629	MIMAT0015642	mmu-miR-3471	
QM043-02	D23	MmiRQP0299	MIMAT0004545	mmu-miR-200b-5p	
QM043-02	D24	MmiRQP1630	MIMAT0015643	mmu-miR-3472	
QM043-02	E01	MmiRQP1387	MIMAT0009455	mmu-miR-1983	
QM043-02	E02	MmiRQP0854	MIMAT0000131	mmu-miR-99a-5p	
QM043-02	E03	MmiRQP1388	MIMAT0009456	mmu-miR-1839-5p	
QM043-02	E04	MmiRQP0862	MIMAT0000385	mmu-miR-106a-5p	
QM043-02	E05	MmiRQP1391	MIMAT0011208	mmu-miR-2132	
QM043-02	E06	MmiRQP0930	MIMAT0000668	mmu-miR-211-5p	
QM043-02	E07	MmiRQP1392	MIMAT0011209	mmu-miR-2133	
QM043-02	E08	MmiRQP0941	MIMAT0000368	mmu-miR-291a-3p	
QM043-02	E09	MmiRQP1393	MIMAT0011210	mmu-miR-2134	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	E10	MmiRQP0951	MIMAT0000375	mmu-miR-297a-5p	
QM043-02	E11	MmiRQP1394	MIMAT0011211	mmu-miR-2135	
QM043-02	E12	MmiRQP0952	MIMAT0003480	mmu-miR-297b-5p	
QM043-02	E13	MmiRQP1395	MIMAT0011212	mmu-miR-2136	
QM043-02	E14	MmiRQP0953	MIMAT0004865	mmu-miR-297c-5p	
QM043-02	E15	MmiRQP1396	MIMAT0011213	mmu-miR-2137	
QM043-02	E16	MmiRQP0959	MIMAT0004186	mmu-miR-301b-3p	
QM043-02	E17	MmiRQP1397	MIMAT0011214	mmu-miR-2138	
QM043-02	E18	MmiRQP1015	MIMAT0004871	mmu-miR-465b-5p	
QM043-02	E19	MmiRQP1398	MIMAT0011215	mmu-miR-2139	
QM043-02	E20	MmiRQP1017	MIMAT0004873	mmu-miR-465c-5p	
QM043-02	E21	MmiRQP1399	MIMAT0011216	mmu-miR-2140	
QM043-02	E22	MmiRQP1018	MIMAT0004759	mmu-miR-466a-5p	
QM043-02	E23	MmiRQP1400	MIMAT0011217	mmu-miR-2141	
QM043-02	E24	MmiRQP1019	MIMAT0005453	mmu-miR-466b-3-3p	
QM043-02	F01	MmiRQP0301	MIMAT0004663	mmu-miR-200c-5p	
QM043-02	F02	MmiRQP1632	MIMAT0015645	mmu-miR-3473	
QM043-02	F03	MmiRQP0328	MIMAT0005444	mmu-miR-218-2-3p	
QM043-02	F04	MmiRQP1633	MIMAT0015646	mmu-miR-3474	
QM043-02	F05	MmiRQP0333	MIMAT0004629	mmu-miR-22-5p	
QM043-02	F06	MmiRQP1643	MIMAT0014816	mmu-miR-3099-3p	
QM043-02	F07	MmiRQP0356	MIMAT0004630	mmu-miR-26b-3p	
QM043-02	F08	MmiRQP1644	MIMAT0014853	mmu-miR-3072-3p	
QM043-02	F09	MmiRQP0358	MIMAT0004633	mmu-miR-27a-5p	
QM043-02	F10	MmiRQP1645	MIMAT0014801	mmu-miR-1247-3p	
QM043-02	F11	MmiRQP0370	MIMAT0004631	mmu-miR-29a-5p	
QM043-02	F12	MmiRQP1646	MIMAT0014802	mmu-miR-1264-5p	
QM043-02	F13	MmiRQP0376	MIMAT0004632	mmu-miR-29c-5p	
QM043-02	F14	MmiRQP1647	MIMAT0014803	mmu-miR-1264-3p	
QM043-02	F15	MmiRQP0390	MIMAT0000129	mmu-miR-30a-3p	
QM043-02	F16	MmiRQP1648	MIMAT0014804	mmu-miR-1249-5p	
QM043-02	F17	MmiRQP0394	MIMAT0004616	mmu-miR-30c-1-3p	
QM043-02	F18	MmiRQP1649	MIMAT0014805	mmu-miR-1843-5p	
QM043-02	F19	MmiRQP0399	MIMAT0000249	mmu-miR-30e-3p	
QM043-02	F20	MmiRQP1650	MIMAT0014806	mmu-miR-1843-3p	
QM043-02	F21	MmiRQP0429	MIMAT0004666	mmu-miR-33-3p	
QM043-02	F22	MmiRQP1651	MIMAT0014807	mmu-miR-344d-3-5p	
QM043-02	F23	MmiRQP0474	MIMAT0000742	mmu-miR-378-5p	
QM043-02	F24	MmiRQP1652	MIMAT0014808	mmu-miR-344d-3p	
QM043-02	G01	MmiRQP1401	MIMAT0011218	mmu-miR-2142	
QM043-02	G02	MmiRQP1022	MIMAT0004931	mmu-miR-466d-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	G03	MmiRQP1402	MIMAT0011219	mmu-miR-2143	
QM043-02	G04	MmiRQP1024	MIMAT0004879	mmu-miR-466e-5p	
QM043-02	G05	MmiRQP1403	MIMAT0011220	mmu-miR-2144	
QM043-02	G06	MmiRQP1027	MIMAT0004881	mmu-miR-466f-5p	
QM043-02	G07	MmiRQP1404	MIMAT0011221	mmu-miR-2145	
QM043-02	G08	MmiRQP1028	MIMAT0005844	mmu-miR-466f	
QM043-02	G09	MmiRQP1405	MIMAT0011222	mmu-miR-2146	
QM043-02	G10	MmiRQP1030	MIMAT0004884	mmu-miR-466h-5p	
QM043-02	G11	MmiRQP1406	MIMAT0011286	mmu-miR-2182	
QM043-02	G12	MmiRQP1032	MIMAT0005848	mmu-miR-466j	
QM043-02	G13	MmiRQP1407	MIMAT0011287	mmu-miR-2183	
QM043-02	G14	MmiRQP1035	MIMAT0003409	mmu-miR-467a-5p	
QM043-02	G15	MmiRQP1410	MIMAT0012773	mmu-miR-767	
QM043-02	G16	MmiRQP1038	MIMAT0004885	mmu-miR-467c-5p	
QM043-02	G17	MmiRQP1411	MIMAT0009397	mmu-miR-1933-3p	
QM043-02	G18	MmiRQP1041	MIMAT0005293	mmu-miR-467e-5p	
QM043-02	G19	MmiRQP1412	MIMAT0009406	mmu-miR-1941-3p	
QM043-02	G20	MmiRQP1045	MIMAT0005855	mmu-miR-467h	
QM043-02	G21	MmiRQP1413	MIMAT0009457	mmu-miR-1839-3p	
QM043-02	G22	MmiRQP1087	MIMAT0005833	mmu-miR-669d-5p	
QM043-02	G23	MmiRQP0007	MIMAT0000383	mmu-let-7d-5p	
QM043-02	G24	MmiRQP1182	MIMAT0004849	mmu-miR-883a-3p	
QM043-02	H01	MmiRQP0516	MIMAT0003485	mmu-miR-455-5p	
QM043-02	H02	MmiRQP1654	MIMAT0014811	mmu-miR-3059-5p	
QM043-02	H03	MmiRQP0784	MIMAT0003498	mmu-miR-708-3p	
QM043-02	H04	MmiRQP1656	MIMAT0014813	mmu-miR-3058-5p	
QM043-02	H05	MmiRQP0785	MIMAT0004670	mmu-miR-7a-1-3p	
QM043-02	H06	MmiRQP1657	MIMAT0014814	mmu-miR-3058-3p	
QM043-02	H07	MmiRQP0789	MIMAT0004820	mmu-miR-744-3p	
QM043-02	H08	MmiRQP1658	MIMAT0014817	mmu-miR-3106-5p	
QM043-02	H09	MmiRQP0824	MIMAT0000143	mmu-miR-9-3p	
QM043-02	H10	MmiRQP1659	MIMAT0014818	mmu-miR-3106-3p	
QM043-02	H11	MmiRQP0836	MIMAT0004636	mmu-miR-93-3p	
QM043-02	H12	MmiRQP1660	MIMAT0014819	mmu-miR-344d-1-5p	
QM043-02	H13	MmiRQP0856	MIMAT0004525	mmu-miR-99b-3p	
QM043-02	H14	MmiRQP1661	MIMAT0014961	mmu-miR-344d-2-5p	
QM043-02	H15	MmiRQP1482	MIMAT0014857	mmu-miR-3074-1-3p	
QM043-02	H16	MmiRQP1662	MIMAT0014822	mmu-miR-3057-5p	
QM043-02	H17	MmiRQP0858	MIMAT0004622	mmu-let-7c-1-3p	
QM043-02	H18	MmiRQP1663	MIMAT0014823	mmu-miR-3057-3p	
QM043-02	H19	MmiRQP0859	MIMAT0005439	mmu-let-7c-2-3p	mmu-let-7a-1-3p

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	H20	MmiRQP1664	MIMAT0014825	mmu-miR-1251-3p	
QM043-02	H21	MmiRQP0860	MIMAT0004519	mmu-let-7g-3p	
QM043-02	H22	MmiRQP1665	MIMAT0014826	mmu-miR-3060-5p	
QM043-02	H23	MmiRQP0863	MIMAT0004538	mmu-miR-10b-3p	
QM043-02	H24	MmiRQP1666	MIMAT0014827	mmu-miR-3060-3p	
QM043-02	I01	MmiRQP0015	MIMAT0000121	mmu-let-7g-5p	
QM043-02	I02	MmiRQP1184	MIMAT0004851	mmu-miR-883b-3p	
QM043-02	I03	MmiRQP0017	MIMAT0000122	mmu-let-7i-5p	
QM043-02	I04	MmiRQP1190	MIMAT0000616	mmu-miR-101b-3p	
QM043-02	I05	MmiRQP0156	MIMAT0000141	mmu-miR-130a-3p	
QM043-02	I06	MmiRQP1287	MIMAT0000678	mmu-miR-7b-5p	
QM043-02	I07	MmiRQP0159	MIMAT0000387	mmu-miR-130b-3p	
QM043-02	I08	MmiRQP1344	MIMAT0009401	mmu-miR-1937a	
QM043-02	I09	MmiRQP0184	MIMAT0000153	mmu-miR-141-3p	
QM043-02	I10	MmiRQP1432	MIMAT0009414	mmu-miR-1937b	
QM043-02	I11	MmiRQP0196	MIMAT0000158	mmu-miR-146a-5p	
QM043-02	I12	MmiRQP1359	MIMAT0009418	mmu-miR-669l-5p	
QM043-02	I13	MmiRQP0197	MIMAT0003475	mmu-miR-146b-5p	
QM043-02	I14	MmiRQP1361	MIMAT0009421	mmu-miR-669o-5p	
QM043-02	I15	MmiRQP0204	MIMAT0000516	mmu-miR-148a-3p	
QM043-02	I16	MmiRQP1369	MIMAT0009429	mmu-miR-1937c	
QM043-02	I17	MmiRQP0206	MIMAT0000580	mmu-miR-148b-3p	
QM043-02	I18	MmiRQP1414	MIMAT0009460	mmu-miR-1982.1-3p	
QM043-02	I19	MmiRQP0255	MIMAT0000528	mmu-miR-18a-5p	
QM043-02	I20	MmiRQP0289	MIMAT0004667	mmu-miR-199b-3p	mmu-miR-199a-3p
QM043-02	I21	MmiRQP0298	MIMAT0000519	mmu-miR-200a-3p	
QM043-02	I22	MmiRQP1367	MIMAT0009427	mmu-miR-669n	
QM043-02	I23	MmiRQP0312	MIMAT0000529	mmu-miR-20a-5p	
QM043-02	I24	MmiRQP1016	MIMAT0004217	mmu-miR-465a-3p	mmu-miR-465c-3p mmu-miR-465b-3p
QM043-02	J01	MmiRQP0879	MIMAT0004583	mmu-miR-130b-5p	
QM043-02	J02	MmiRQP1667	MIMAT0014828	mmu-miR-3061-5p	
QM043-02	J03	MmiRQP0883	MIMAT0004668	mmu-miR-138-1-3p	
QM043-02	J04	MmiRQP1668	MIMAT0014829	mmu-miR-3061-3p	
QM043-02	J05	MmiRQP0885	MIMAT0004533	mmu-miR-141-5p	
QM043-02	J06	MmiRQP1669	MIMAT0014830	mmu-miR-3062-5p	
QM043-02	J07	MmiRQP0886	MIMAT0004534	mmu-miR-145-3p	
QM043-02	J08	MmiRQP1670	MIMAT0014831	mmu-miR-3062-3p	
QM043-02	J09	MmiRQP0887	MIMAT0004826	mmu-miR-146b-3p	
QM043-02	J10	MmiRQP1671	MIMAT0014832	mmu-miR-3063-5p	
QM043-02	J11	MmiRQP0891	MIMAT0004624	mmu-miR-15a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	J12	MmiRQP1672	MIMAT0014833	mmu-miR-3063-3p	
QM043-02	J13	MmiRQP0892	MIMAT0004625	mmu-miR-16-1-3p	
QM043-02	J14	MmiRQP1673	MIMAT0014834	mmu-miR-3064-5p	
QM043-02	J15	MmiRQP0893	MIMAT0000650	mmu-miR-17-3p	
QM043-02	J16	MmiRQP1676	MIMAT0014837	mmu-miR-3065-3p	
QM043-02	J17	MmiRQP0894	MIMAT0005443	mmu-miR-181a-2-3p	
QM043-02	J18	MmiRQP1677	MIMAT0014838	mmu-miR-3066-5p	
QM043-02	J19	MmiRQP0896	MIMAT0004540	mmu-miR-186-3p	
QM043-02	J20	MmiRQP1678	MIMAT0014839	mmu-miR-3066-3p	
QM043-02	J21	MmiRQP0907	MIMAT0004626	mmu-miR-18a-3p	
QM043-02	J22	MmiRQP1679	MIMAT0014840	mmu-miR-3067-5p	
QM043-02	J23	MmiRQP0917	MIMAT0004542	mmu-miR-191-3p	
QM043-02	J24	MmiRQP1681	MIMAT0014842	mmu-miR-3068-5p	
QM043-02	K01	MmiRQP0314	MIMAT0003187	mmu-miR-20b-5p	
QM043-02	K02	MmiRQP1025	MIMAT0002107	mmu-miR-466a-3p	mmu-miR-466e-3p mmu-miR-466c-3p mmu-miR-466b-3p
QM043-02	K03	MmiRQP0387	MIMAT0003377	mmu-miR-302d-3p	
QM043-02	K04	MmiRQP0110	MIMAT0004530	mmu-miR-127-5p	
QM043-02	K05	MmiRQP0444	MIMAT0000381	mmu-miR-34c-5p	
QM043-02	K06	MmiRQP0882	MIMAT0004532	mmu-miR-136-3p	
QM043-02	K07	MmiRQP0835	MIMAT0004899	mmu-miR-92b-3p	
QM043-02	K08	MmiRQP0218	MIMAT0004537	mmu-miR-154-3p	
QM043-02	K09	MmiRQP0853	MIMAT0000545	mmu-miR-98-5p	
QM043-02	K10	MmiRQP0369	MIMAT0000377	mmu-miR-299-5p	
QM043-02	K11	MmiRQP0908	MIMAT0004858	mmu-miR-18b-5p	
QM043-02	K12	MmiRQP1223	MIMAT0004578	mmu-miR-300-5p	
QM043-02	K13	MmiRQP0980	MIMAT0000382	mmu-miR-34b-5p	
QM043-02	K14	MmiRQP0988	MIMAT0003387	mmu-miR-376a-5p	
QM043-02	K15	MmiRQP1040	MIMAT0004886	mmu-miR-467d-5p	
QM043-02	K16	MmiRQP1254	MIMAT0003388	mmu-miR-376b-5p	
QM043-02	K17	MmiRQP1044	MIMAT0005854	mmu-miR-467g	
QM043-02	K18	MmiRQP0991	MIMAT0005295	mmu-miR-376c-5p	
QM043-02	K19	MmiRQP1085	MIMAT0003476	mmu-miR-669b-5p	
QM043-02	K20	MmiRQP0994	MIMAT0004691	mmu-miR-382-3p	
QM043-02	K21	MmiRQP1086	MIMAT0003479	mmu-miR-669c-5p	
QM043-02	K22	MmiRQP0487	MIMAT0001093	mmu-miR-411-3p	
QM043-02	K23	MmiRQP1092	MIMAT0005842	mmu-miR-669h-3p	
QM043-02	K24	MmiRQP0498	MIMAT0004753	mmu-miR-431-3p	
QM043-02	L01	MmiRQP0919	MIMAT0004618	mmu-miR-196a-2-3p	
QM043-02	L02	MmiRQP1682	MIMAT0014843	mmu-miR-3068-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	L03	MmiRQP0920	MIMAT0000672	mmu-miR-199b-5p	
QM043-02	L04	MmiRQP1684	MIMAT0014845	mmu-miR-3069-3p	
QM043-02	L05	MmiRQP0921	MIMAT0004660	mmu-miR-19a-5p	
QM043-02	L06	MmiRQP1685	MIMAT0014846	mmu-miR-3070a-5p	mmu-miR-3070b-5p
QM043-02	L07	MmiRQP0925	MIMAT0004547	mmu-miR-203-5p	
QM043-02	L08	MmiRQP1686	MIMAT0014847	mmu-miR-3070a-3p	
QM043-02	L09	MmiRQP0927	MIMAT0004627	mmu-miR-20a-3p	
QM043-02	L10	MmiRQP1687	MIMAT0014849	mmu-miR-3070b-3p	
QM043-02	L11	MmiRQP0934	MIMAT0004665	mmu-miR-218-1-3p	
QM043-02	L12	MmiRQP1688	MIMAT0014850	mmu-miR-3071-5p	
QM043-02	L13	MmiRQP0937	MIMAT0000218	mmu-miR-24-1-5p	
QM043-02	L14	MmiRQP1690	MIMAT0014854	mmu-miR-3073-5p	
QM043-02	L15	MmiRQP0938	MIMAT0004661	mmu-miR-28-3p	
QM043-02	L16	MmiRQP1694	MIMAT0014859	mmu-miR-3075-3p	
QM043-02	L17	MmiRQP0945	MIMAT0004573	mmu-miR-293-5p	
QM043-02	L18	MmiRQP1695	MIMAT0014860	mmu-miR-3076-5p	
QM043-02	L19	MmiRQP0948	MIMAT0004574	mmu-miR-294-5p	
QM043-02	L20	MmiRQP1696	MIMAT0014861	mmu-miR-3076-3p	
QM043-02	L21	MmiRQP0950	MIMAT0004575	mmu-miR-295-5p	
QM043-02	L22	MmiRQP1698	MIMAT0014863	mmu-miR-3077-3p	
QM043-02	L23	MmiRQP0957	MIMAT0004523	mmu-miR-29b-1-5p	
QM043-02	L24	MmiRQP1699	MIMAT0014864	mmu-miR-3078-5p	
QM043-02	M01	MmiRQP1293	MIMAT0000539	mmu-miR-92a-3p	
QM043-02	M02	MmiRQP1000	MIMAT0001419	mmu-miR-433-5p	
QM043-02	M03	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM043-02	M04	MmiRQP1053	MIMAT0003129	mmu-miR-485-3p	
QM043-02	M05	MmiRQP0004	MIMAT0000522	mmu-let-7b-5p	
QM043-02	M06	MmiRQP1627	MIMAT0015640	mmu-miR-3470a	
QM043-02	M07	MmiRQP0006	MIMAT0000523	mmu-let-7c-5p	
QM043-02	M08	MmiRQP1631	MIMAT0015644	mmu-miR-1186b	
QM043-02	M09	MmiRQP0010	MIMAT0000524	mmu-let-7e-5p	
QM043-02	M10	MmiRQP0877	MIMAT0004529	mmu-miR-125b-2-3p	
QM043-02	M11	MmiRQP0012	MIMAT0000525	mmu-let-7f-5p	
QM043-02	M12	MmiRQP2749	MIMAT0017282	mmu-miR-544-5p	
QM043-02	M13	MmiRQP0018	MIMAT0000655	mmu-miR-100-5p	
QM043-02	M14	MmiRQP0880	MIMAT0003473	mmu-miR-133a-5p	
QM043-02	M15	MmiRQP0021	MIMAT0000133	mmu-miR-101a-3p	
QM043-02	M16	MmiRQP2750	MIMAT0017254	mmu-miR-499-3p	
QM043-02	M17	MmiRQP0022	MIMAT0000546	mmu-miR-103-3p	
QM043-02	M18	MmiRQP2219	MIMAT0020367	mmu-miR-3473b	
QM043-02	M19	MmiRQP0030	MIMAT0000647	mmu-miR-107-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	M20	MmiRQP1062	MIMAT0004790	mmu-miR-503-3p	
QM043-02	M21	MmiRQP0032	MIMAT0000648	mmu-miR-10a-5p	
QM043-02	M22	MmiRQP2568	MIMAT0017022	mmu-miR-34a-3p	
QM043-02	M23	MmiRQP0034	MIMAT0000208	mmu-miR-10b-5p	
QM043-02	M24	MmiRQP0929	MIMAT0004628	mmu-miR-21-3p	
QM043-02	N01	MmiRQP0960	MIMAT0004579	mmu-miR-302a-5p	
QM043-02	N02	MmiRQP1700	MIMAT0014865	mmu-miR-3078-3p	
QM043-02	N03	MmiRQP0961	MIMAT0003373	mmu-miR-302b-5p	
QM043-02	N04	MmiRQP1702	MIMAT0014867	mmu-miR-3079-3p	
QM043-02	N05	MmiRQP0963	MIMAT0003375	mmu-miR-302c-5p	
QM043-02	N06	MmiRQP1704	MIMAT0014869	mmu-miR-3080-3p	
QM043-02	N07	MmiRQP0964	MIMAT0004524	mmu-miR-30b-3p	
QM043-02	N08	MmiRQP1705	MIMAT0014870	mmu-miR-3081-5p	
QM043-02	N09	MmiRQP0965	MIMAT0004634	mmu-miR-31-3p	
QM043-02	N10	MmiRQP1706	MIMAT0014871	mmu-miR-3081-3p	
QM043-02	N11	MmiRQP0966	MIMAT0000549	mmu-miR-322-3p	
QM043-02	N12	MmiRQP1707	MIMAT0014872	mmu-miR-3082-5p	
QM043-02	N13	MmiRQP0970	MIMAT0000569	mmu-miR-330-3p	
QM043-02	N14	MmiRQP1708	MIMAT0014873	mmu-miR-3082-3p	
QM043-02	N15	MmiRQP0981	MIMAT0004580	mmu-miR-34c-3p	
QM043-02	N16	MmiRQP1710	MIMAT0014875	mmu-miR-3083-3p	
QM043-02	N17	MmiRQP0986	MIMAT0003728	mmu-miR-374-3p	
QM043-02	N18	MmiRQP1712	MIMAT0014877	mmu-miR-3084-3p	
QM043-02	N19	MmiRQP0998	MIMAT0001342	mmu-miR-425-3p	
QM043-02	N20	MmiRQP1713	MIMAT0014878	mmu-miR-3085-5p	
QM043-02	N21	MmiRQP1011	MIMAT0002104	mmu-miR-463-5p	
QM043-02	N22	MmiRQP1714	MIMAT0014879	mmu-miR-3085-3p	
QM043-02	N23	MmiRQP1036	MIMAT0003478	mmu-miR-467b-3p	
QM043-02	N24	MmiRQP1715	MIMAT0014880	mmu-miR-3086-5p	
QM043-02	O01	MmiRQP2685	MIMAT0022365	mmu-miR-5619-5p	
QM043-02	O02	MmiRQP2698	MIMAT0022378	mmu-miR-5624-3p	
QM043-02	O03	MmiRQP2686	MIMAT0022366	mmu-miR-5619-3p	
QM043-02	O04	MmiRQP2699	MIMAT0022379	mmu-miR-5625-5p	
QM043-02	O05	MmiRQP2687	MIMAT0022367	mmu-miR-5620-5p	
QM043-02	O06	MmiRQP2700	MIMAT0022380	mmu-miR-5625-3p	
QM043-02	O07	MmiRQP2688	MIMAT0022368	mmu-miR-5620-3p	
QM043-02	O08	MmiRQP2702	MIMAT0022382	mmu-miR-5626-3p	
QM043-02	O09	MmiRQP2689	MIMAT0022369	mmu-miR-5621-5p	
QM043-02	O10	MmiRQP2703	MIMAT0022385	mmu-miR-5627-5p	
QM043-02	O11	MmiRQP2691	MIMAT0022371	mmu-miR-5622-5p	
QM043-02	O12	MmiRQP2704	MIMAT0022386	mmu-miR-5627-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-02	O13	MmiRQP2692	MIMAT0022372	mmu-miR-5622-3p	
QM043-02	O14	MmiRQP2740	MIMAT0022503	mmu-miR-344i	
QM043-02	O15	MmiRQP2693	MIMAT0022373	mmu-miR-5623-5p	
QM043-02	O16	MmiRQP2741	MIMAT0022504	mmu-miR-5709	
QM043-02	O17	MmiRQP2694	MIMAT0022374	mmu-miR-5623-3p	
QM043-02	O18	MmiRQP2742	MIMAT0022505	mmu-miR-5710	
QM043-02	O19	MmiRQP2695	MIMAT0022375	mmu-miR-3073b-5p	
QM043-02	O20	MmiRQP2746	MIMAT0017345	mmu-miR-669l-3p	
QM043-02	O21	MmiRQP2696	MIMAT0022376	mmu-miR-3073b-3p	
QM043-02	O22	MmiRQP2747	MIMAT0017047	mmu-miR-1a-2-5p	
QM043-02	O23	MmiRQP2697	MIMAT0022377	mmu-miR-5624-5p	
QM043-02	O24	MmiRQP2748	MIMAT0017324	mmu-miR-669d-3p	
QM043-02	P01	NC	NC	NC	
QM043-02	P02	NC	NC	NC	
QM043-02	P03	NC	NC	NC	
QM043-02	P04	NC	NC	NC	
QM043-02	P05	MmiRQP9002	NR_003027	MK1	
QM043-02	P06	MmiRQP9002	NR_003027	MK1	
QM043-02	P07	MmiRQP9012	NR_028128	MK2	
QM043-02	P08	MmiRQP9012	NR_028128	MK2	
QM043-02	P09	MmiRQP9022	NR_028554	MK3	
QM043-02	P10	MmiRQP9022	NR_028554	MK3	
QM043-02	P11	MmiRQP9042	NR_028550	MK4	
QM043-02	P12	MmiRQP9042	NR_028550	MK4	
QM043-02	P13	MmiRQP9052	NR_028541	MK5	
QM043-02	P14	MmiRQP9052	NR_028541	MK5	
QM043-02	P15	MmiRQP9062	NR_028091	MK6	
QM043-02	P16	MmiRQP9062	NR_028091	MK6	
QM043-02	P17	RT	RT	RT	
QM043-02	P18	RT	RT	RT	
QM043-02	P19	RT	RT	RT	
QM043-02	P20	RT	RT	RT	
QM043-02	P21	PCR	PCR	PCR	
QM043-02	P22	PCR	PCR	PCR	
QM043-02	P23	PCR	PCR	PCR	
QM043-02	P24	PCR	PCR	PCR	
QM043-03	A01	MmiRQP1716	MIMAT0014881	mmu-miR-3086-3p	
QM043-03	A02	MmiRQP2534	MIMAT0020614	mmu-miR-3473c	
QM043-03	A03	MmiRQP1717	MIMAT0014882	mmu-miR-466m-5p	
QM043-03	A04	MmiRQP2535	MIMAT0020615	mmu-miR-5107	
QM043-03	A05	MmiRQP1718	MIMAT0014883	mmu-miR-466m-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	A06	MmiRQP2536	MIMAT0020616	mmu-miR-5108	
QM043-03	A07	MmiRQP1719	MIMAT0014884	mmu-miR-669d-2-3p	
QM043-03	A08	MmiRQP2537	MIMAT0020617	mmu-miR-5109	
QM043-03	A09	MmiRQP1720	MIMAT0014885	mmu-miR-466o-5p	mmu-miR-466b-5p
QM043-03	A10	MmiRQP2538	MIMAT0020618	mmu-miR-5110	
QM043-03	A11	MmiRQP1721	MIMAT0014886	mmu-miR-466o-3p	
QM043-03	A12	MmiRQP2539	MIMAT0020619	mmu-miR-5111-3p	
QM043-03	A13	MmiRQP1722	MIMAT0002108	mmu-miR-467a-3p	mmu-miR-467d-3p
QM043-03	A14	MmiRQP2540	MIMAT0020620	mmu-miR-5112	
QM043-03	A15	MmiRQP1724	MIMAT0017243	mmu-miR-669a-3p	
QM043-03	A16	MmiRQP2541	MIMAT0020621	mmu-miR-5113	
QM043-03	A17	MmiRQP1727	MIMAT0014890	mmu-miR-669p-3p	
QM043-03	A18	MmiRQP2542	MIMAT0020622	mmu-miR-5114	
QM043-03	A19	MmiRQP1728	MIMAT0014891	mmu-miR-466p-5p	
QM043-03	A20	MmiRQP2543	MIMAT0020623	mmu-miR-5115	
QM043-03	A21	MmiRQP1729	MIMAT0014893	mmu-miR-466n-5p	
QM043-03	A22	MmiRQP2544	MIMAT0020624	mmu-miR-5116	
QM043-03	A23	MmiRQP1730	MIMAT0014894	mmu-miR-466n-3p	
QM043-03	A24	MmiRQP2545	MIMAT0020625	mmu-miR-5117-5p	
QM043-03	B01				
QM043-03	B02				
QM043-03	B03				
QM043-03	B04				
QM043-03	B05				
QM043-03	B06				
QM043-03	B07				
QM043-03	B08				
QM043-03	B09				
QM043-03	B10				
QM043-03	B11				
QM043-03	B12				
QM043-03	B13				
QM043-03	B14				
QM043-03	B15				
QM043-03	B16				
QM043-03	B17				
QM043-03	B18				
QM043-03	B19				
QM043-03	B20				
QM043-03	B21				
QM043-03	B22				

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	B23				
QM043-03	B24				
QM043-03	C01	MmiRQP1731	MIMAT0014895	mmu-miR-3087-5p	
QM043-03	C02	MmiRQP2546	MIMAT0020626	mmu-miR-5118	
QM043-03	C03	MmiRQP1732	MIMAT0014896	mmu-miR-3087-3p	
QM043-03	C04	MmiRQP2550	MIMAT0020630	mmu-miR-5122	
QM043-03	C05	MmiRQP1733	MIMAT0014897	mmu-miR-3088-5p	
QM043-03	C06	MmiRQP2551	MIMAT0020631	mmu-miR-466q	
QM043-03	C07	MmiRQP1734	MIMAT0014898	mmu-miR-3088-3p	
QM043-03	C08	MmiRQP2552	MIMAT0020632	mmu-miR-3473d	
QM043-03	C09	MmiRQP1735	MIMAT0014899	mmu-miR-3089-5p	
QM043-03	C10	MmiRQP2554	MIMAT0020634	mmu-miR-5124	
QM043-03	C11	MmiRQP1736	MIMAT0014900	mmu-miR-3089-3p	
QM043-03	C12	MmiRQP2555	MIMAT0020635	mmu-miR-5125	
QM043-03	C13	MmiRQP1737	MIMAT0014901	mmu-miR-3090-5p	
QM043-03	C14				
QM043-03	C15	MmiRQP1738	MIMAT0014902	mmu-miR-3090-3p	
QM043-03	C16				
QM043-03	C17	MmiRQP1739	MIMAT0014903	mmu-miR-3091-5p	
QM043-03	C18				
QM043-03	C19	MmiRQP1740	MIMAT0014904	mmu-miR-3091-3p	
QM043-03	C20				
QM043-03	C21	MmiRQP1741	MIMAT0014905	mmu-miR-3092-5p	
QM043-03	C22				
QM043-03	C23	MmiRQP1742	MIMAT0014906	mmu-miR-3092-3p	
QM043-03	C24				
QM043-03	D01				
QM043-03	D02				
QM043-03	D03				
QM043-03	D04				
QM043-03	D05				
QM043-03	D06				
QM043-03	D07				
QM043-03	D08				
QM043-03	D09				
QM043-03	D10				
QM043-03	D11				
QM043-03	D12				
QM043-03	D13				
QM043-03	D14				
QM043-03	D15				

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	D16				
QM043-03	D17				
QM043-03	D18				
QM043-03	D19				
QM043-03	D20				
QM043-03	D21				
QM043-03	D22				
QM043-03	D23				
QM043-03	D24				
QM043-03	E01	MmiRQP1743	MIMAT0014907	mmu-miR-3093-5p	
QM043-03	E02				
QM043-03	E03	MmiRQP1745	MIMAT0014909	mmu-miR-3094-5p	
QM043-03	E04				
QM043-03	E05	MmiRQP1746	MIMAT0014910	mmu-miR-3094-3p	
QM043-03	E06				
QM043-03	E07	MmiRQP1747	MIMAT0014911	mmu-miR-3095-5p	
QM043-03	E08				
QM043-03	E09	MmiRQP1748	MIMAT0014912	mmu-miR-3095-3p	
QM043-03	E10				
QM043-03	E11	MmiRQP1749	MIMAT0014913	mmu-miR-3096-5p	
QM043-03	E12				
QM043-03	E13	MmiRQP1750	MIMAT0014914	mmu-miR-3096-3p	
QM043-03	E14				
QM043-03	E15	MmiRQP1751	MIMAT0014915	mmu-miR-3097-5p	
QM043-03	E16				
QM043-03	E17	MmiRQP1753	MIMAT0014917	mmu-miR-3098-5p	
QM043-03	E18				
QM043-03	E19	MmiRQP1756	MIMAT0014920	mmu-miR-3100-3p	
QM043-03	E20				
QM043-03	E21	MmiRQP1758	MIMAT0014922	mmu-miR-3101-3p	
QM043-03	E22				
QM043-03	E23	MmiRQP1759	MIMAT0014923	mmu-miR-344e-5p	mmu-miR-344h-5p
QM043-03	E24				
QM043-03	F01				
QM043-03	F02				
QM043-03	F03				
QM043-03	F04				
QM043-03	F05				
QM043-03	F06				
QM043-03	F07				
QM043-03	F08				

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	F09				
QM043-03	F10				
QM043-03	F11				
QM043-03	F12				
QM043-03	F13				
QM043-03	F14				
QM043-03	F15				
QM043-03	F16				
QM043-03	F17				
QM043-03	F18				
QM043-03	F19				
QM043-03	F20				
QM043-03	F21				
QM043-03	F22				
QM043-03	F23				
QM043-03	F24				
QM043-03	G01	MmiRQP1761	MIMAT0014925	mmu-miR-344b-5p	
QM043-03	G02				
QM043-03	G03	MmiRQP1762	MIMAT0014926	mmu-miR-344b-3p	
QM043-03	G04				
QM043-03	G05	MmiRQP1764	MIMAT0014928	mmu-miR-344c-3p	
QM043-03	G06				
QM043-03	G07	MmiRQP1765	MIMAT0014929	mmu-miR-344g-5p	
QM043-03	G08				
QM043-03	G09	MmiRQP1766	MIMAT0014930	mmu-miR-344g-3p	
QM043-03	G10				
QM043-03	G11	MmiRQP1767	MIMAT0014931	mmu-miR-344f-5p	
QM043-03	G12				
QM043-03	G13	MmiRQP1768	MIMAT0014932	mmu-miR-344f-3p	
QM043-03	G14				
QM043-03	G15	MmiRQP1770	MIMAT0014934	mmu-miR-3102-5p.2-5p	
QM043-03	G16				
QM043-03	G17	MmiRQP1771	MIMAT0014935	mmu-miR-3102-3p.2	
QM043-03	G18				
QM043-03	G19	MmiRQP1772	MIMAT0014936	mmu-miR-3102	
QM043-03	G20				
QM043-03	G21	MmiRQP1773	MIMAT0014937	mmu-miR-3103-5p	
QM043-03	G22				
QM043-03	G23	MmiRQP1774	MIMAT0014938	mmu-miR-3103-3p	
QM043-03	G24				
QM043-03	H01				

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	H02				
QM043-03	H03				
QM043-03	H04				
QM043-03	H05				
QM043-03	H06				
QM043-03	H07				
QM043-03	H08				
QM043-03	H09				
QM043-03	H10				
QM043-03	H11				
QM043-03	H12				
QM043-03	H13				
QM043-03	H14				
QM043-03	H15				
QM043-03	H16				
QM043-03	H17				
QM043-03	H18				
QM043-03	H19				
QM043-03	H20				
QM043-03	H21				
QM043-03	H22				
QM043-03	H23				
QM043-03	H24				
QM043-03	I01	MmiRQP1775	MIMAT0014939	mmu-miR-3104-5p	
QM043-03	I02				
QM043-03	I03	MmiRQP1776	MIMAT0014940	mmu-miR-3104-3p	
QM043-03	I04				
QM043-03	I05	MmiRQP1777	MIMAT0014941	mmu-miR-3105-5p	
QM043-03	I06				
QM043-03	I07	MmiRQP1778	MIMAT0014942	mmu-miR-3105-3p	
QM043-03	I08				
QM043-03	I09	MmiRQP1779	MIMAT0014944	mmu-miR-3107-3p	
QM043-03	I10				
QM043-03	I11	MmiRQP1780	MIMAT0014946	mmu-miR-3074-2-3p	
QM043-03	I12				
QM043-03	I13	MmiRQP1781	MIMAT0014947	mmu-miR-3108-5p	
QM043-03	I14				
QM043-03	I15	MmiRQP1782	MIMAT0014948	mmu-miR-3108-3p	
QM043-03	I16				
QM043-03	I17	MmiRQP1783	MIMAT0014949	mmu-miR-3109-5p	
QM043-03	I18				

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	I19	MmiRQP1784	MIMAT0014950	mmu-miR-3109-3p	
QM043-03	I20				
QM043-03	I21	MmiRQP1785	MIMAT0014951	mmu-miR-3110-5p	
QM043-03	I22				
QM043-03	I23	MmiRQP1786	MIMAT0014952	mmu-miR-3110-3p	
QM043-03	I24				
QM043-03	J01				
QM043-03	J02				
QM043-03	J03				
QM043-03	J04				
QM043-03	J05				
QM043-03	J06				
QM043-03	J07				
QM043-03	J08				
QM043-03	J09				
QM043-03	J10				
QM043-03	J11				
QM043-03	J12				
QM043-03	J13				
QM043-03	J14				
QM043-03	J15				
QM043-03	J16				
QM043-03	J17				
QM043-03	J18				
QM043-03	J19				
QM043-03	J20				
QM043-03	J21				
QM043-03	J22				
QM043-03	J23				
QM043-03	J24				
QM043-03	K01	MmiRQP1787	MIMAT0014953	mmu-miR-374c-5p	
QM043-03	K02				
QM043-03	K03	MmiRQP1788	MIMAT0014954	mmu-miR-374c-3p	
QM043-03	K04				
QM043-03	K05	MmiRQP1789	MIMAT0014955	mmu-miR-3112-5p	
QM043-03	K06				
QM043-03	K07	MmiRQP1790	MIMAT0014956	mmu-miR-3112-3p	
QM043-03	K08				
QM043-03	K09	MmiRQP1791	MIMAT0014957	mmu-miR-1912-5p	
QM043-03	K10				
QM043-03	K11	MmiRQP1793	MIMAT0014959	mmu-miR-3113-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	K12				
QM043-03	K13	MmiRQP1810	MIMAT0022384	mmu-miR-344h-3p	
QM043-03	K14				
QM043-03	K15	MmiRQP2192	MIMAT0019336	mmu-miR-3960	
QM043-03	K16				
QM043-03	K17	MmiRQP2193	MIMAT0019338	mmu-miR-3961	
QM043-03	K18				
QM043-03	K19	MmiRQP2195	MIMAT0019340	mmu-miR-3962	
QM043-03	K20				
QM043-03	K21	MmiRQP2196	MIMAT0019341	mmu-miR-3963	
QM043-03	K22				
QM043-03	K23	MmiRQP2197	MIMAT0019342	mmu-miR-3096b-5p	
QM043-03	K24				
QM043-03	L01				
QM043-03	L02				
QM043-03	L03				
QM043-03	L04				
QM043-03	L05				
QM043-03	L06				
QM043-03	L07				
QM043-03	L08				
QM043-03	L09				
QM043-03	L10				
QM043-03	L11				
QM043-03	L12				
QM043-03	L13				
QM043-03	L14				
QM043-03	L15				
QM043-03	L16				
QM043-03	L17				
QM043-03	L18				
QM043-03	L19				
QM043-03	L20				
QM043-03	L21				
QM043-03	L22				
QM043-03	L23				
QM043-03	L24				
QM043-03	M01	MmiRQP2198	MIMAT0019343	mmu-miR-3096b-3p	
QM043-03	M02				
QM043-03	M03	MmiRQP2199	MIMAT0019344	mmu-miR-3964	
QM043-03	M04				

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	M05	MmiRQP2201	MIMAT0019346	mmu-miR-1843b-3p	
QM043-03	M06				
QM043-03	M07	MmiRQP2202	MIMAT0019347	mmu-miR-3965	
QM043-03	M08				
QM043-03	M09	MmiRQP2203	MIMAT0019348	mmu-miR-378b	
QM043-03	M10				
QM043-03	M11	MmiRQP2204	MIMAT0019349	mmu-miR-101c	
QM043-03	M12				
QM043-03	M13	MmiRQP2205	MIMAT0019350	mmu-miR-3966	
QM043-03	M14				
QM043-03	M15	MmiRQP2206	MIMAT0019351	mmu-miR-3967	
QM043-03	M16				
QM043-03	M17	MmiRQP2207	MIMAT0019352	mmu-miR-3968	
QM043-03	M18				
QM043-03	M19	MmiRQP2208	MIMAT0019353	mmu-miR-3969	
QM043-03	M20				
QM043-03	M21	MmiRQP2209	MIMAT0019354	mmu-miR-28b	
QM043-03	M22				
QM043-03	M23	MmiRQP2210	MIMAT0019355	mmu-miR-3970	
QM043-03	M24				
QM043-03	N01				
QM043-03	N02				
QM043-03	N03				
QM043-03	N04				
QM043-03	N05				
QM043-03	N06				
QM043-03	N07				
QM043-03	N08				
QM043-03	N09				
QM043-03	N10				
QM043-03	N11				
QM043-03	N12				
QM043-03	N13				
QM043-03	N14				
QM043-03	N15				
QM043-03	N16				
QM043-03	N17				
QM043-03	N18				
QM043-03	N19				
QM043-03	N20				
QM043-03	N21				

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	N22				
QM043-03	N23				
QM043-03	N24				
QM043-03	O01	MmiRQP2211	MIMAT0019356	mmu-miR-3971	
QM043-03	O02				
QM043-03	O03	MmiRQP2518	MIMAT0020540	mmu-miR-5046	
QM043-03	O04				
QM043-03	O05	MmiRQP2524	MIMAT0020604	mmu-miR-5097	
QM043-03	O06				
QM043-03	O07	MmiRQP2525	MIMAT0020605	mmu-miR-5098	
QM043-03	O08				
QM043-03	O09	MmiRQP2526	MIMAT0020606	mmu-miR-5099	
QM043-03	O10				
QM043-03	O11	MmiRQP2527	MIMAT0020607	mmu-miR-5100	
QM043-03	O12				
QM043-03	O13	MmiRQP2528	MIMAT0020608	mmu-miR-5101	
QM043-03	O14				
QM043-03	O15	MmiRQP2529	MIMAT0020609	mmu-miR-5102	
QM043-03	O16				
QM043-03	O17	MmiRQP2530	MIMAT0020610	mmu-miR-5103	
QM043-03	O18				
QM043-03	O19	MmiRQP2531	MIMAT0020611	mmu-miR-5104	
QM043-03	O20				
QM043-03	O21	MmiRQP2532	MIMAT0020612	mmu-miR-5105	
QM043-03	O22				
QM043-03	O23	MmiRQP2533	MIMAT0020613	mmu-miR-5106	
QM043-03	O24				
QM043-03	P01	NC	NC	NC	
QM043-03	P02	NC	NC	NC	
QM043-03	P03	NC	NC	NC	
QM043-03	P04	NC	NC	NC	
QM043-03	P05	MmiRQP9002	NR_003027	MK1	
QM043-03	P06	MmiRQP9002	NR_003027	MK1	
QM043-03	P07	MmiRQP9012	NR_028128	MK2	
QM043-03	P08	MmiRQP9012	NR_028128	MK2	
QM043-03	P09	MmiRQP9022	NR_028554	MK3	
QM043-03	P10	MmiRQP9022	NR_028554	MK3	
QM043-03	P11	MmiRQP9042	NR_028550	MK4	
QM043-03	P12	MmiRQP9042	NR_028550	MK4	
QM043-03	P13	MmiRQP9052	NR_028541	MK5	
QM043-03	P14	MmiRQP9052	NR_028541	MK5	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM043-03	P15	MmiRQP9062	NR_028091	MK6	
QM043-03	P16	MmiRQP9062	NR_028091	MK6	
QM043-03	P17	RT	RT	RT	
QM043-03	P18	RT	RT	RT	
QM043-03	P19	RT	RT	RT	
QM043-03	P20	RT	RT	RT	
QM043-03	P21	PCR	PCR	PCR	
QM043-03	P22	PCR	PCR	PCR	
QM043-03	P23	PCR	PCR	PCR	
QM043-03	P24	PCR	PCR	PCR	