

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	A01	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM045_01	A02	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM045_01	A03	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM045_01	A04	HmiRQP0006	MIMAT0000064	hsa-let-7c-5p	
QM045_01	A05	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM045_01	A06	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM045_01	A07	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM045_01	A08	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM045_01	A09	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM045_01	A10	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM045_01	A11	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM045_01	A12	HmiRQP0019	MIMAT0004512	hsa-miR-100-3p	
QM045_01	A13	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM045_01	A14	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM045_01	A15	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM045_01	A16	HmiRQP4680	MIMAT0037306	hsa-miR-103a-1-5p	
QM045_01	A17	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM045_01	A18	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM045_01	A19	HmiRQP0023	MIMAT0007402	hsa-miR-103b	
QM045_01	A20	HmiRQP0025	MIMAT0000102	hsa-miR-105-5p	
QM045_01	A21	HmiRQP0027	MIMAT0004517	hsa-miR-106a-3p	
QM045_01	A22	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM045_01	A23	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM045_01	A24	HmiRQP0029	MIMAT0000680	hsa-miR-106b-5p	
QM045_01	B01	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM045_01	B02	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM045_01	B03	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM045_01	B04	HmiRQP0033	MIMAT0004556	hsa-miR-10b-3p	
QM045_01	B05	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM045_01	B06	HmiRQP3286	MIMAT0022713	hsa-miR-1185-2-3p	
QM045_01	B07	HmiRQP0054	MIMAT0005873	hsa-miR-1208	
QM045_01	B08	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM045_01	B09	HmiRQP0057	MIMAT0005459	hsa-miR-1224-3p	
QM045_01	B10	HmiRQP0058	MIMAT0005458	hsa-miR-1224-5p	
QM045_01	B11	HmiRQP0060	MIMAT0005573	hsa-miR-1225-3p	
QM045_01	B12	HmiRQP0059	MIMAT0005572	hsa-miR-1225-5p	
QM045_01	B13	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM045_01	B14	HmiRQP0062	MIMAT0005577	hsa-miR-1226-3p	
QM045_01	B15	HmiRQP0061	MIMAT0005576	hsa-miR-1226-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	B16	HmiRQP0063	MIMAT0005580	hsa-miR-1227-3p	
QM045_01	B17	HmiRQP3394	MIMAT0022941	hsa-miR-1227-5p	
QM045_01	B18	HmiRQP0065	MIMAT0005583	hsa-miR-1228-3p	
QM045_01	B19	HmiRQP0064	MIMAT0005582	hsa-miR-1228-5p	
QM045_01	B20	HmiRQP0066	MIMAT0005584	hsa-miR-1229-3p	
QM045_01	B21	HmiRQP0070	MIMAT0005591	hsa-miR-1236-3p	
QM045_01	B22	HmiRQP3397	MIMAT0022945	hsa-miR-1236-5p	
QM045_01	B23	HmiRQP0072	MIMAT0005593	hsa-miR-1238-3p	
QM045_01	B24	HmiRQP3399	MIMAT0022947	hsa-miR-1238-5p	
QM045_01	C01	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM045_01	C02	HmiRQP0078	MIMAT0005898	hsa-miR-1246	
QM045_01	C03	HmiRQP3407	MIMAT0022721	hsa-miR-1247-3p	
QM045_01	C04	HmiRQP0087	MIMAT0005906	hsa-miR-1255a	
QM045_01	C05	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM045_01	C06	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM045_01	C07	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM045_01	C08	HmiRQP0097	MIMAT0004603	hsa-miR-125b-2-3p	
QM045_01	C09	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM045_01	C10	HmiRQP0100	MIMAT0005911	hsa-miR-1260a	
QM045_01	C11	HmiRQP1496	MIMAT0015041	hsa-miR-1260b	
QM045_01	C12	HmiRQP0101	MIMAT0005913	hsa-miR-1261	
QM045_01	C13	HmiRQP0102	MIMAT0005914	hsa-miR-1262	
QM045_01	C14	HmiRQP0103	MIMAT0005915	hsa-miR-1263	
QM045_01	C15	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM045_01	C16	HmiRQP0104	MIMAT0005791	hsa-miR-1264	
QM045_01	C17	HmiRQP0105	MIMAT0005918	hsa-miR-1265	
QM045_01	C18	HmiRQP0098	MIMAT0000444	hsa-miR-126-5p	
QM045_01	C19	HmiRQP3410	MIMAT0026742	hsa-miR-1266-3p	
QM045_01	C20	HmiRQP0106	MIMAT0005920	hsa-miR-1266-5p	
QM045_01	C21	HmiRQP0107	MIMAT0005921	hsa-miR-1267	
QM045_01	C22	HmiRQP0108	MIMAT0005922	hsa-miR-1268a	
QM045_01	C23	HmiRQP2034	MIMAT0018925	hsa-miR-1268b	
QM045_01	C24	HmiRQP0109	MIMAT0005923	hsa-miR-1269a	
QM045_01	D01	HmiRQP2166	MIMAT0019059	hsa-miR-1269b	
QM045_01	D02	HmiRQP3284	MIMAT0022712	hsa-miR-1271-3p	
QM045_01	D03	HmiRQP0113	MIMAT0005796	hsa-miR-1271-5p	
QM045_01	D04	HmiRQP0115	MIMAT0005926	hsa-miR-1273a	
QM045_01	D05	HmiRQP2521	MIMAT0020601	hsa-miR-1273f	
QM045_01	D06	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	D07	HmiRQP0118	MIMAT0005929	hsa-miR-1275	
QM045_01	D08	HmiRQP0125	MIMAT0000424	hsa-miR-128-3p	
QM045_01	D09	HmiRQP0134	MIMAT0005880	hsa-miR-1290	
QM045_01	D10	HmiRQP0139	MIMAT0004605	hsa-miR-129-2-3p	
QM045_01	D11	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM045_01	D12	HmiRQP0152	MIMAT0005893	hsa-miR-1305	
QM045_01	D13	HmiRQP0154	MIMAT0005951	hsa-miR-1307-3p	
QM045_01	D14	HmiRQP3416	MIMAT0022727	hsa-miR-1307-5p	
QM045_01	D15	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM045_01	D16	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM045_01	D17	HmiRQP0159	MIMAT0000691	hsa-miR-130b-3p	
QM045_01	D18	HmiRQP0158	MIMAT0004680	hsa-miR-130b-5p	
QM045_01	D19	HmiRQP0163	MIMAT0005953	hsa-miR-1322	
QM045_01	D20	HmiRQP0164	MIMAT0005795	hsa-miR-1323	
QM045_01	D21	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM045_01	D22	HmiRQP0165	MIMAT0005956	hsa-miR-1324	
QM045_01	D23	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM045_01	D24	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM045_01	E01	HmiRQP2307	MIMAT0019776	hsa-miR-1343-3p	
QM045_01	E02	HmiRQP3060	MIMAT0026481	hsa-miR-134-3p	
QM045_01	E03	HmiRQP0168	MIMAT0000447	hsa-miR-134-5p	
QM045_01	E04	HmiRQP0170	MIMAT0000428	hsa-miR-135a-5p	
QM045_01	E05	HmiRQP0171	MIMAT0004698	hsa-miR-135b-3p	
QM045_01	E06	HmiRQP0172	MIMAT0000758	hsa-miR-135b-5p	
QM045_01	E07	HmiRQP0174	MIMAT0004606	hsa-miR-136-3p	
QM045_01	E08	HmiRQP0173	MIMAT0000448	hsa-miR-136-5p	
QM045_01	E09	HmiRQP0175	MIMAT0000429	hsa-miR-137	
QM045_01	E10	HmiRQP4720	MIMAT0037310	hsa-miR-137-5p	
QM045_01	E11	HmiRQP0177	MIMAT0000430	hsa-miR-138-5p	
QM045_01	E12	HmiRQP3046	MIMAT0000250	hsa-miR-139-5p	
QM045_01	E13	HmiRQP0044	MIMAT0000416	hsa-miR-1-3p	
QM045_01	E14	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM045_01	E15	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM045_01	E16	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM045_01	E17	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM045_01	E18	HmiRQP0186	MIMAT0000434	hsa-miR-142-3p	
QM045_01	E19	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM045_01	E20	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM045_01	E21	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	E22	HmiRQP0190	MIMAT0000436	hsa-miR-144-3p	
QM045_01	E23	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM045_01	E24	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM045_01	F01	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM045_01	F02	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM045_01	F03	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM045_01	F04	HmiRQP0198	MIMAT0004766	hsa-miR-146b-3p	
QM045_01	F05	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM045_01	F06	HmiRQP0204	MIMAT0000243	hsa-miR-148a-3p	
QM045_01	F07	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM045_01	F08	HmiRQP0206	MIMAT0000759	hsa-miR-148b-3p	
QM045_01	F09	HmiRQP0205	MIMAT0004699	hsa-miR-148b-5p	
QM045_01	F10	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM045_01	F11	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM045_01	F12	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM045_01	F13	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM045_01	F14	HmiRQP0211	MIMAT0000757	hsa-miR-151a-3p	
QM045_01	F15	HmiRQP0212	MIMAT0004697	hsa-miR-151a-5p	
QM045_01	F16	HmiRQP0219	MIMAT0000452	hsa-miR-154-5p	
QM045_01	F17	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM045_01	F18	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM045_01	F19	HmiRQP2185	MIMAT0019077	hsa-miR-1587	
QM045_01	F20	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM045_01	F21	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM045_01	F22	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM045_01	F23	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM045_01	F24	HmiRQP0228	MIMAT0004518	hsa-miR-16-2-3p	
QM045_01	G01	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM045_01	G02	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM045_01	G03	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM045_01	G04	HmiRQP0233	MIMAT0004558	hsa-miR-181a-2-3p	
QM045_01	G05	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM045_01	G06	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM045_01	G07	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM045_01	G08	HmiRQP0236	MIMAT0004559	hsa-miR-181c-3p	
QM045_01	G09	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM045_01	G10	HmiRQP0237	MIMAT0002821	hsa-miR-181d-5p	
QM045_01	G11	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM045_01	G12	HmiRQP0240	MIMAT0006765	hsa-miR-1825	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	G13	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM045_01	G14	HmiRQP0242	MIMAT0006767	hsa-miR-1827	
QM045_01	G15	HmiRQP0243	MIMAT0004560	hsa-miR-183-3p	
QM045_01	G16	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM045_01	G17	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM045_01	G18	HmiRQP0246	MIMAT0004611	hsa-miR-185-3p	
QM045_01	G19	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM045_01	G20	HmiRQP0249	MIMAT0004612	hsa-miR-186-3p	
QM045_01	G21	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM045_01	G22	HmiRQP0251	MIMAT0000262	hsa-miR-187-3p	
QM045_01	G23	HmiRQP0254	MIMAT0002891	hsa-miR-18a-3p	
QM045_01	G24	HmiRQP0255	MIMAT0000072	hsa-miR-18a-5p	
QM045_01	H01	HmiRQP0265	MIMAT0007884	hsa-miR-1910-5p	
QM045_01	H02	HmiRQP0266	MIMAT0007886	hsa-miR-1911-3p	
QM045_01	H03	HmiRQP0268	MIMAT0007887	hsa-miR-1912	
QM045_01	H04	HmiRQP0269	MIMAT0007888	hsa-miR-1913	
QM045_01	H05	HmiRQP0264	MIMAT0001618	hsa-miR-191-3p	
QM045_01	H06	HmiRQP0270	MIMAT0007889	hsa-miR-1914-5p	
QM045_01	H07	HmiRQP0273	MIMAT0007892	hsa-miR-1915-3p	
QM045_01	H08	HmiRQP0272	MIMAT0007891	hsa-miR-1915-5p	
QM045_01	H09	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM045_01	H10	HmiRQP0275	MIMAT0004543	hsa-miR-192-3p	
QM045_01	H11	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM045_01	H12	HmiRQP0276	MIMAT0000459	hsa-miR-193a-3p	
QM045_01	H13	HmiRQP0277	MIMAT0004614	hsa-miR-193a-5p	
QM045_01	H14	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM045_01	H15	HmiRQP0279	MIMAT0004767	hsa-miR-193b-5p	
QM045_01	H16	HmiRQP0280	MIMAT0000460	hsa-miR-194-5p	
QM045_01	H17	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM045_01	H18	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM045_01	H19	HmiRQP0285	MIMAT0004562	hsa-miR-196a-3p	
QM045_01	H20	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM045_01	H21	HmiRQP0286	MIMAT0001080	hsa-miR-196b-5p	
QM045_01	H22	HmiRQP0287	MIMAT0000227	hsa-miR-197-3p	
QM045_01	H23	HmiRQP3038	MIMAT0022691	hsa-miR-197-5p	
QM045_01	H24	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM045_01	I01	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM045_01	I02	HmiRQP0289	MIMAT0004563	hsa-miR-199b-3p	
QM045_01	I03	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	I04	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM045_01	I05	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM045_01	I06	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM045_01	I07	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	
QM045_01	I08	HmiRQP0297	MIMAT0001620	hsa-miR-200a-5p	
QM045_01	I09	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM045_01	I10	HmiRQP0299	MIMAT0004571	hsa-miR-200b-5p	
QM045_01	I11	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM045_01	I12	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM045_01	I13	HmiRQP0305	MIMAT0000264	hsa-miR-203a-3p	
QM045_01	I14	HmiRQP2344	MIMAT0019814	hsa-miR-203b-3p	
QM045_01	I15	HmiRQP2343	MIMAT0019813	hsa-miR-203b-5p	
QM045_01	I16	HmiRQP3040	MIMAT0022693	hsa-miR-204-3p	
QM045_01	I17	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM045_01	I18	HmiRQP1331	MIMAT0009977	hsa-miR-2052	
QM045_01	I19	HmiRQP1314	MIMAT0009197	hsa-miR-205-3p	
QM045_01	I20	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM045_01	I21	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM045_01	I22	HmiRQP0309	MIMAT0000241	hsa-miR-208a-3p	
QM045_01	I23	HmiRQP3043	MIMAT0026474	hsa-miR-208a-5p	
QM045_01	I24	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM045_01	J01	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM045_01	J02	HmiRQP0313	MIMAT0004752	hsa-miR-20b-3p	
QM045_01	J03	HmiRQP0314	MIMAT0001413	hsa-miR-20b-5p	
QM045_01	J04	HmiRQP0317	MIMAT0000267	hsa-miR-210-3p	
QM045_01	J05	HmiRQP3049	MIMAT0026475	hsa-miR-210-5p	
QM045_01	J06	HmiRQP1333	MIMAT0010133	hsa-miR-2110	
QM045_01	J07	HmiRQP3016	MIMAT0022694	hsa-miR-211-3p	
QM045_01	J08	HmiRQP1306	MIMAT0011157	hsa-miR-2114-3p	
QM045_01	J09	HmiRQP1324	MIMAT0011156	hsa-miR-2114-5p	
QM045_01	J10	HmiRQP1302	MIMAT0011159	hsa-miR-2115-3p	
QM045_01	J11	HmiRQP1297	MIMAT0011158	hsa-miR-2115-5p	
QM045_01	J12	HmiRQP0318	MIMAT0000268	hsa-miR-211-5p	
QM045_01	J13	HmiRQP1304	MIMAT0011161	hsa-miR-2116-3p	
QM045_01	J14	HmiRQP1319	MIMAT0011160	hsa-miR-2116-5p	
QM045_01	J15	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM045_01	J16	HmiRQP3050	MIMAT0022695	hsa-miR-212-5p	
QM045_01	J17	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM045_01	J18	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	



Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	J19	HmiRQP0321	MIMAT0004564	hsa-miR-214-5p	
QM045_01	J20	HmiRQP3051	MIMAT0026476	hsa-miR-215-3p	
QM045_01	J21	HmiRQP0322	MIMAT0000272	hsa-miR-215-5p	
QM045_01	J22	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM045_01	J23	HmiRQP3052	MIMAT0022844	hsa-miR-216a-3p	
QM045_01	J24	HmiRQP0323	MIMAT0000273	hsa-miR-216a-5p	
QM045_01	K01	HmiRQP0324	MIMAT0004959	hsa-miR-216b-5p	
QM045_01	K02	HmiRQP0325	MIMAT0000274	hsa-miR-217	
QM045_01	K03	HmiRQP0326	MIMAT0004565	hsa-miR-218-1-3p	
QM045_01	K04	HmiRQP0328	MIMAT0004566	hsa-miR-218-2-3p	
QM045_01	K05	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM045_01	K06	HmiRQP0329	MIMAT0004567	hsa-miR-219a-1-3p	
QM045_01	K07	HmiRQP0330	MIMAT0000276	hsa-miR-219a-5p	
QM045_01	K08	HmiRQP1876	MIMAT0019747	hsa-miR-219b-5p	
QM045_01	K09	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM045_01	K10	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM045_01	K11	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM045_01	K12	HmiRQP0340	MIMAT0004569	hsa-miR-222-5p	
QM045_01	K13	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM045_01	K14	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM045_01	K15	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM045_01	K16	HmiRQP1295	MIMAT0009198	hsa-miR-224-3p	
QM045_01	K17	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM045_01	K18	HmiRQP0333	MIMAT0004495	hsa-miR-22-5p	
QM045_01	K19	HmiRQP1327	MIMAT0011775	hsa-miR-2276-3p	
QM045_01	K20	HmiRQP1328	MIMAT0011777	hsa-miR-2277-3p	
QM045_01	K21	HmiRQP3434	MIMAT0017352	hsa-miR-2277-5p	
QM045_01	K22	HmiRQP1312	MIMAT0011778	hsa-miR-2278	
QM045_01	K23	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM045_01	K24	HmiRQP0345	MIMAT0004496	hsa-miR-23a-5p	
QM045_01	L01	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM045_01	L02	HmiRQP0347	MIMAT0004587	hsa-miR-23b-5p	
QM045_01	L03	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM045_01	L04	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM045_01	L05	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM045_01	L06	HmiRQP0355	MIMAT0004681	hsa-miR-26a-2-3p	
QM045_01	L07	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM045_01	L08	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM045_01	L09	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	L10	HmiRQP0358	MIMAT0004501	hsa-miR-27a-5p	
QM045_01	L11	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM045_01	L12	HmiRQP0360	MIMAT0004588	hsa-miR-27b-5p	
QM045_01	L13	HmiRQP0363	MIMAT0004502	hsa-miR-28-3p	
QM045_01	L14	HmiRQP0362	MIMAT0000085	hsa-miR-28-5p	
QM045_01	L15	HmiRQP0364	MIMAT0000690	hsa-miR-296-5p	
QM045_01	L16	HmiRQP0368	MIMAT0000687	hsa-miR-299-3p	
QM045_01	L17	HmiRQP0369	MIMAT0002890	hsa-miR-299-5p	
QM045_01	L18	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM045_01	L19	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM045_01	L20	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM045_01	L21	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM045_01	L22	HmiRQP0376	MIMAT0004673	hsa-miR-29c-5p	
QM045_01	L23	HmiRQP0377	MIMAT0004903	hsa-miR-300	
QM045_01	L24	HmiRQP0378	MIMAT0000688	hsa-miR-301a-3p	
QM045_01	M01	HmiRQP0383	MIMAT0000715	hsa-miR-302b-3p	
QM045_01	M02	HmiRQP1518	MIMAT0015066	hsa-miR-3065-5p	
QM045_01	M03	HmiRQP0390	MIMAT0000088	hsa-miR-30a-3p	
QM045_01	M04	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM045_01	M05	HmiRQP0392	MIMAT0004589	hsa-miR-30b-3p	
QM045_01	M06	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM045_01	M07	HmiRQP0394	MIMAT0004674	hsa-miR-30c-1-3p	
QM045_01	M08	HmiRQP0395	MIMAT0004550	hsa-miR-30c-2-3p	
QM045_01	M09	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM045_01	M10	HmiRQP0397	MIMAT0004551	hsa-miR-30d-3p	
QM045_01	M11	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM045_01	M12	HmiRQP0399	MIMAT0000693	hsa-miR-30e-3p	
QM045_01	M13	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM045_01	M14	HmiRQP1435	MIMAT0014977	hsa-miR-3115	
QM045_01	M15	HmiRQP1436	MIMAT0014978	hsa-miR-3116	
QM045_01	M16	HmiRQP1437	MIMAT0014979	hsa-miR-3117-3p	
QM045_01	M17	HmiRQP1439	MIMAT0014981	hsa-miR-3119	
QM045_01	M18	HmiRQP1440	MIMAT0014982	hsa-miR-3120-3p	
QM045_01	M19	HmiRQP1441	MIMAT0014983	hsa-miR-3121-3p	
QM045_01	M20	HmiRQP1442	MIMAT0014984	hsa-miR-3122	
QM045_01	M21	HmiRQP1444	MIMAT0014986	hsa-miR-3124-5p	
QM045_01	M22	HmiRQP1446	MIMAT0014988	hsa-miR-3125	
QM045_01	M23	HmiRQP1634	MIMAT0015377	hsa-miR-3126-3p	
QM045_01	M24	HmiRQP1447	MIMAT0014989	hsa-miR-3126-5p	



Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	N01	HmiRQP1448	MIMAT0014990	hsa-miR-3127-5p	
QM045_01	N02	HmiRQP1450	MIMAT0014992	hsa-miR-3129-5p	
QM045_01	N03	HmiRQP1635	MIMAT0014994	hsa-miR-3130-3p	
QM045_01	N04	HmiRQP1451	MIMAT0014995	hsa-miR-3130-5p	
QM045_01	N05	HmiRQP1452	MIMAT0014996	hsa-miR-3131	
QM045_01	N06	HmiRQP1453	MIMAT0014997	hsa-miR-3132	
QM045_01	N07	HmiRQP1454	MIMAT0014998	hsa-miR-3133	
QM045_01	N08	HmiRQP1456	MIMAT0015000	hsa-miR-3134	
QM045_01	N09	HmiRQP1457	MIMAT0015001	hsa-miR-3135a	
QM045_01	N10	HmiRQP2093	MIMAT0018985	hsa-miR-3135b	
QM045_01	N11	HmiRQP1459	MIMAT0015003	hsa-miR-3136-5p	
QM045_01	N12	HmiRQP1461	MIMAT0015005	hsa-miR-3137	
QM045_01	N13	HmiRQP1462	MIMAT0015006	hsa-miR-3138	
QM045_01	N14	HmiRQP1463	MIMAT0015007	hsa-miR-3139	
QM045_01	N15	HmiRQP0402	MIMAT0004504	hsa-miR-31-3p	
QM045_01	N16	HmiRQP1464	MIMAT0015008	hsa-miR-3140-3p	
QM045_01	N17	HmiRQP1466	MIMAT0015010	hsa-miR-3141	
QM045_01	N18	HmiRQP1468	MIMAT0015012	hsa-miR-3143	
QM045_01	N19	HmiRQP1636	MIMAT0015015	hsa-miR-3144-3p	
QM045_01	N20	HmiRQP1471	MIMAT0015016	hsa-miR-3145-3p	
QM045_01	N21	HmiRQP1473	MIMAT0015018	hsa-miR-3146	
QM045_01	N22	HmiRQP1474	MIMAT0015019	hsa-miR-3147	
QM045_01	N23	HmiRQP1476	MIMAT0015021	hsa-miR-3148	
QM045_01	N24	HmiRQP1477	MIMAT0015022	hsa-miR-3149	
QM045_01	O01	HmiRQP1478	MIMAT0015023	hsa-miR-3150a-3p	
QM045_01	O02	HmiRQP3475	MIMAT0027026	hsa-miR-3151-3p	
QM045_01	O03	HmiRQP1479	MIMAT0015024	hsa-miR-3151-5p	
QM045_01	O04	HmiRQP1480	MIMAT0015025	hsa-miR-3152-3p	
QM045_01	O05	HmiRQP1481	MIMAT0015026	hsa-miR-3153	
QM045_01	O06	HmiRQP1483	MIMAT0015028	hsa-miR-3154	
QM045_01	O07	HmiRQP1484	MIMAT0015029	hsa-miR-3155a	
QM045_01	O08	HmiRQP3013	MIMAT0019209	hsa-miR-3156-3p	
QM045_01	O09	HmiRQP1485	MIMAT0015030	hsa-miR-3156-5p	
QM045_01	O10	HmiRQP1486	MIMAT0015031	hsa-miR-3157-5p	
QM045_01	O11	HmiRQP1487	MIMAT0015032	hsa-miR-3158-3p	
QM045_01	O12	HmiRQP1488	MIMAT0015033	hsa-miR-3159	
QM045_01	O13	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM045_01	O14	HmiRQP1489	MIMAT0015034	hsa-miR-3160-3p	
QM045_01	O15	HmiRQP1490	MIMAT0015035	hsa-miR-3161	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_01	O16	HmiRQP3481	MIMAT0019213	hsa-miR-3162-3p	
QM045_01	O17	HmiRQP1491	MIMAT0015036	hsa-miR-3162-5p	
QM045_01	O18	HmiRQP1492	MIMAT0015037	hsa-miR-3163	
QM045_01	O19	HmiRQP1493	MIMAT0015038	hsa-miR-3164	
QM045_01	O20	HmiRQP1494	MIMAT0015039	hsa-miR-3165	
QM045_01	O21	HmiRQP1495	MIMAT0015040	hsa-miR-3166	
QM045_01	O22	HmiRQP1497	MIMAT0015042	hsa-miR-3167	
QM045_01	O23	HmiRQP1498	MIMAT0015043	hsa-miR-3168	
QM045_01	O24	HmiRQP1499	MIMAT0015044	hsa-miR-3169	
QM045_01	P01	NC	NC	NC	
QM045_01	P02	NC	NC	NC	
QM045_01	P03	NC	NC	NC	
QM045_01	P04	NC	NC	NC	
QM045_01	P05	HmiRQP9001	NR_002752	HK1	
QM045_01	P06	HmiRQP9001	NR_002752	HK1	
QM045_01	P07	HmiRQP9011	NR_002750	HK2	
QM045_01	P08	HmiRQP9011	NR_002750	HK2	
QM045_01	P09	HmiRQP9021	NR_002745	HK3	
QM045_01	P10	HmiRQP9021	NR_002745	HK3	
QM045_01	P11	HmiRQP9051	NR_002746	HK4	
QM045_01	P12	HmiRQP9051	NR_002746	HK4	
QM045_01	P13	HmiRQP9061	NR_002744	HK5	
QM045_01	P14	HmiRQP9061	NR_002744	HK5	
QM045_01	P15	HmiRQP9071	NR_002450	HK6	
QM045_01	P16	HmiRQP9071	NR_002450	HK6	
QM045_01	P17	RT	RT	RT	
QM045_01	P18	RT	RT	RT	
QM045_01	P19	RT	RT	RT	
QM045_01	P20	RT	RT	RT	
QM045_01	P21	PCR	PCR	PCR	
QM045_01	P22	PCR	PCR	PCR	
QM045_01	P23	PCR	PCR	PCR	
QM045_01	P24	PCR	PCR	PCR	
QM045_02	A01	HmiRQP1500	MIMAT0015045	hsa-miR-3170	
QM045_02	A02	HmiRQP1503	MIMAT0015048	hsa-miR-3173-3p	
QM045_02	A03	HmiRQP1505	MIMAT0015051	hsa-miR-3174	
QM045_02	A04	HmiRQP1506	MIMAT0015052	hsa-miR-3175	
QM045_02	A05	HmiRQP1507	MIMAT0015053	hsa-miR-3176	
QM045_02	A06	HmiRQP1508	MIMAT0015054	hsa-miR-3177-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_02	A07	HmiRQP1509	MIMAT0015055	hsa-miR-3178	
QM045_02	A08	HmiRQP1510	MIMAT0015056	hsa-miR-3179	
QM045_02	A09	HmiRQP1982	MIMAT0018178	hsa-miR-3180	
QM045_02	A10	HmiRQP1638	MIMAT0015058	hsa-miR-3180-3p	
QM045_02	A11	HmiRQP1511	MIMAT0015057	hsa-miR-3180-5p	
QM045_02	A12	HmiRQP1513	MIMAT0015061	hsa-miR-3181	
QM045_02	A13	HmiRQP1514	MIMAT0015062	hsa-miR-3182	
QM045_02	A14	HmiRQP1515	MIMAT0015063	hsa-miR-3183	
QM045_02	A15	HmiRQP3484	MIMAT0022731	hsa-miR-3184-3p	
QM045_02	A16	HmiRQP1516	MIMAT0015064	hsa-miR-3184-5p	
QM045_02	A17	HmiRQP1517	MIMAT0015065	hsa-miR-3185	
QM045_02	A18	HmiRQP1640	MIMAT0015068	hsa-miR-3186-3p	
QM045_02	A19	HmiRQP1519	MIMAT0015067	hsa-miR-3186-5p	
QM045_02	A20	HmiRQP1520	MIMAT0015069	hsa-miR-3187-3p	
QM045_02	A21	HmiRQP3485	MIMAT0019216	hsa-miR-3187-5p	
QM045_02	A22	HmiRQP1521	MIMAT0015070	hsa-miR-3188	
QM045_02	A23	HmiRQP1522	MIMAT0015071	hsa-miR-3189-3p	
QM045_02	A24	HmiRQP3487	MIMAT0015073	hsa-miR-3190-5p	
QM045_02	B01	HmiRQP1525	MIMAT0015075	hsa-miR-3191-3p	
QM045_02	B02	HmiRQP3489	MIMAT0027027	hsa-miR-3192-3p	
QM045_02	B03	HmiRQP1526	MIMAT0015076	hsa-miR-3192-5p	
QM045_02	B04	HmiRQP1527	MIMAT0015077	hsa-miR-3193	
QM045_02	B05	HmiRQP3490	MIMAT0019218	hsa-miR-3194-3p	
QM045_02	B06	HmiRQP1528	MIMAT0015078	hsa-miR-3194-5p	
QM045_02	B07	HmiRQP1529	MIMAT0015079	hsa-miR-3195	
QM045_02	B08	HmiRQP1530	MIMAT0015080	hsa-miR-3196	
QM045_02	B09	HmiRQP1532	MIMAT0015082	hsa-miR-3197	
QM045_02	B10	HmiRQP1533	MIMAT0015083	hsa-miR-3198	
QM045_02	B11	HmiRQP1534	MIMAT0015084	hsa-miR-3199	
QM045_02	B12	HmiRQP0405	MIMAT0000510	hsa-miR-320a	
QM045_02	B13	HmiRQP4731	MIMAT0037311	hsa-miR-320a-5p	
QM045_02	B14	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM045_02	B15	HmiRQP0407	MIMAT0005793	hsa-miR-320c	
QM045_02	B16	HmiRQP1523	MIMAT0015072	hsa-miR-320e	
QM045_02	B17	HmiRQP0410	MIMAT0000755	hsa-miR-323a-3p	
QM045_02	B18	HmiRQP1637	MIMAT0015050	hsa-miR-323b-3p	
QM045_02	B19	HmiRQP0404	MIMAT0000090	hsa-miR-32-5p	
QM045_02	B20	HmiRQP0414	MIMAT0000756	hsa-miR-326	
QM045_02	B21	HmiRQP0415	MIMAT0000752	hsa-miR-328-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_02	B22	HmiRQP0417	MIMAT0000751	hsa-miR-330-3p	
QM045_02	B23	HmiRQP0418	MIMAT0004693	hsa-miR-330-5p	
QM045_02	B24	HmiRQP0420	MIMAT0000760	hsa-miR-331-3p	
QM045_02	C01	HmiRQP0422	MIMAT0004703	hsa-miR-335-3p	
QM045_02	C02	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM045_02	C03	HmiRQP0423	MIMAT0000754	hsa-miR-337-3p	
QM045_02	C04	HmiRQP0424	MIMAT0004695	hsa-miR-337-5p	
QM045_02	C05	HmiRQP0426	MIMAT0000763	hsa-miR-338-3p	
QM045_02	C06	HmiRQP0425	MIMAT0004701	hsa-miR-338-5p	
QM045_02	C07	HmiRQP0428	MIMAT0004702	hsa-miR-339-3p	
QM045_02	C08	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM045_02	C09	HmiRQP0434	MIMAT0004692	hsa-miR-340-5p	
QM045_02	C10	HmiRQP0436	MIMAT0000753	hsa-miR-342-3p	
QM045_02	C11	HmiRQP0435	MIMAT0004694	hsa-miR-342-5p	
QM045_02	C12	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM045_02	C13	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM045_02	C14	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM045_02	C15	HmiRQP0442	MIMAT0000685	hsa-miR-34b-5p	
QM045_02	C16	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM045_02	C17	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	
QM045_02	C18	HmiRQP2408	MIMAT0019877	hsa-miR-3591-3p	
QM045_02	C19	HmiRQP2407	MIMAT0019876	hsa-miR-3591-5p	
QM045_02	C20	HmiRQP1907	MIMAT0017991	hsa-miR-3613-3p	
QM045_02	C21	HmiRQP1910	MIMAT0017994	hsa-miR-3615	
QM045_02	C22	HmiRQP0446	MIMAT0000703	hsa-miR-361-5p	
QM045_02	C23	HmiRQP0448	MIMAT0000705	hsa-miR-362-5p	
QM045_02	C24	HmiRQP0449	MIMAT0000707	hsa-miR-363-3p	
QM045_02	D01	HmiRQP0450	MIMAT0003385	hsa-miR-363-5p	
QM045_02	D02	HmiRQP0451	MIMAT0000710	hsa-miR-365a-3p	
QM045_02	D03	HmiRQP1944	MIMAT0018087	hsa-miR-3665	
QM045_02	D04	HmiRQP0454	MIMAT0000721	hsa-miR-369-3p	
QM045_02	D05	HmiRQP0456	MIMAT0000722	hsa-miR-370-3p	
QM045_02	D06	HmiRQP0458	MIMAT0004687	hsa-miR-371a-5p	
QM045_02	D07	HmiRQP0463	MIMAT0000727	hsa-miR-374a-5p	
QM045_02	D08	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM045_02	D09	HmiRQP0466	MIMAT0000728	hsa-miR-375	
QM045_02	D10	HmiRQP4732	MIMAT0037313	hsa-miR-375-5p	
QM045_02	D11	HmiRQP0467	MIMAT0000729	hsa-miR-376a-3p	
QM045_02	D12	HmiRQP0472	MIMAT0000730	hsa-miR-377-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_02	D13	HmiRQP3114	MIMAT0000732	hsa-miR-378a-3p	
QM045_02	D14	HmiRQP2046	MIMAT0018937	hsa-miR-378g	
QM045_02	D15	HmiRQP2182	MIMAT0019074	hsa-miR-378i	
QM045_02	D16	HmiRQP0475	MIMAT0004690	hsa-miR-379-3p	
QM045_02	D17	HmiRQP0476	MIMAT0000733	hsa-miR-379-5p	
QM045_02	D18	HmiRQP0480	MIMAT0000737	hsa-miR-382-5p	
QM045_02	D19	HmiRQP2192	MIMAT0019337	hsa-miR-3960	
QM045_02	D20	HmiRQP0484	MIMAT0001639	hsa-miR-409-3p	
QM045_02	D21	HmiRQP0483	MIMAT0001638	hsa-miR-409-5p	
QM045_02	D22	HmiRQP0485	MIMAT0002171	hsa-miR-410-3p	
QM045_02	D23	HmiRQP0487	MIMAT0004813	hsa-miR-411-3p	
QM045_02	D24	HmiRQP0486	MIMAT0003329	hsa-miR-411-5p	
QM045_02	E01	HmiRQP0490	MIMAT0001339	hsa-miR-422a	
QM045_02	E02	HmiRQP0491	MIMAT0001340	hsa-miR-423-3p	
QM045_02	E03	HmiRQP0492	MIMAT0004748	hsa-miR-423-5p	
QM045_02	E04	HmiRQP0493	MIMAT0004749	hsa-miR-424-3p	
QM045_02	E05	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM045_02	E06	HmiRQP0496	MIMAT0001343	hsa-miR-425-3p	
QM045_02	E07	HmiRQP0495	MIMAT0003393	hsa-miR-425-5p	
QM045_02	E08	HmiRQP1574	MIMAT0016878	hsa-miR-4257	
QM045_02	E09	HmiRQP1611	MIMAT0016915	hsa-miR-4284	
QM045_02	E10	HmiRQP1612	MIMAT0016916	hsa-miR-4286	
QM045_02	E11	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM045_02	E12	HmiRQP1554	MIMAT0016858	hsa-miR-4306	
QM045_02	E13	HmiRQP0499	MIMAT0001625	hsa-miR-431-5p	
QM045_02	E14	HmiRQP1569	MIMAT0016873	hsa-miR-4322	
QM045_02	E15	HmiRQP0501	MIMAT0002814	hsa-miR-432-5p	
QM045_02	E16	HmiRQP2076	MIMAT0018967	hsa-miR-4448	
QM045_02	E17	HmiRQP2084	MIMAT0018976	hsa-miR-4454	
QM045_02	E18	HmiRQP2100	MIMAT0018993	hsa-miR-4466	
QM045_02	E19	HmiRQP0503	MIMAT0001532	hsa-miR-448	
QM045_02	E20	HmiRQP2124	MIMAT0019018	hsa-miR-4484	
QM045_02	E21	HmiRQP2128	MIMAT0019022	hsa-miR-4488	
QM045_02	E22	HmiRQP0504	MIMAT0001541	hsa-miR-449a	
QM045_02	E23	HmiRQP2147	MIMAT0019041	hsa-miR-4505	
QM045_02	E24	HmiRQP0506	MIMAT0001545	hsa-miR-450a-5p	
QM045_02	F01	HmiRQP0508	MIMAT0004909	hsa-miR-450b-5p	
QM045_02	F02	HmiRQP2159	MIMAT0019053	hsa-miR-4516	
QM045_02	F03	HmiRQP0509	MIMAT0001631	hsa-miR-451a	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_02	F04	HmiRQP2171	MIMAT0019064	hsa-miR-4525	
QM045_02	F05	HmiRQP0510	MIMAT0001635	hsa-miR-452-5p	
QM045_02	F06	HmiRQP0514	MIMAT0003885	hsa-miR-454-3p	
QM045_02	F07	HmiRQP0516	MIMAT0003150	hsa-miR-455-5p	
QM045_02	F08	HmiRQP2231	MIMAT0019699	hsa-miR-4640-5p	
QM045_02	F09	HmiRQP2236	MIMAT0019704	hsa-miR-4644	
QM045_02	F10	HmiRQP2241	MIMAT0019709	hsa-miR-4647	
QM045_02	F11	HmiRQP2275	MIMAT0019743	hsa-miR-4667-5p	
QM045_02	F12	HmiRQP2277	MIMAT0019745	hsa-miR-4668-5p	
QM045_02	F13	HmiRQP2402	MIMAT0019871	hsa-miR-4741	
QM045_02	F14	HmiRQP2431	MIMAT0019900	hsa-miR-4756-3p	
QM045_02	F15	HmiRQP2430	MIMAT0019899	hsa-miR-4756-5p	
QM045_02	F16	HmiRQP2450	MIMAT0019919	hsa-miR-4767	
QM045_02	F17	HmiRQP2458	MIMAT0019927	hsa-miR-4772-3p	
QM045_02	F18	HmiRQP2495	MIMAT0019964	hsa-miR-4792	
QM045_02	F19	HmiRQP0518	MIMAT0002173	hsa-miR-483-3p	
QM045_02	F20	HmiRQP0517	MIMAT0004761	hsa-miR-483-5p	
QM045_02	F21	HmiRQP0519	MIMAT0002174	hsa-miR-484	
QM045_02	F22	HmiRQP0523	MIMAT0002177	hsa-miR-486-5p	
QM045_02	F23	HmiRQP0535	MIMAT0002813	hsa-miR-493-5p	
QM045_02	F24	HmiRQP0536	MIMAT0002816	hsa-miR-494-3p	
QM045_02	G01	HmiRQP0537	MIMAT0002817	hsa-miR-495-3p	
QM045_02	G02	HmiRQP3211	MIMAT0022924	hsa-miR-495-5p	
QM045_02	G03	HmiRQP0540	MIMAT0002820	hsa-miR-497-5p	
QM045_02	G04	HmiRQP0543	MIMAT0002870	hsa-miR-499a-5p	
QM045_02	G05	HmiRQP0544	MIMAT0002871	hsa-miR-500a-3p	
QM045_02	G06	HmiRQP2595	MIMAT0021044	hsa-miR-5010-3p	
QM045_02	G07	HmiRQP2594	MIMAT0021043	hsa-miR-5010-5p	
QM045_02	G08	HmiRQP2597	MIMAT0021046	hsa-miR-5011-3p	
QM045_02	G09	HmiRQP2596	MIMAT0021045	hsa-miR-5011-5p	
QM045_02	G10	HmiRQP0547	MIMAT0004774	hsa-miR-501-3p	
QM045_02	G11	HmiRQP0546	MIMAT0002872	hsa-miR-501-5p	
QM045_02	G12	HmiRQP0548	MIMAT0004775	hsa-miR-502-3p	
QM045_02	G13	HmiRQP0550	MIMAT0002874	hsa-miR-503-5p	
QM045_02	G14	HmiRQP2523	MIMAT0020603	hsa-miR-5096	
QM045_02	G15	HmiRQP2626	MIMAT0022259	hsa-miR-5100	
QM045_02	G16	HmiRQP1537	MIMAT0015087	hsa-miR-514b-5p	
QM045_02	G17	HmiRQP0610	MIMAT0002838	hsa-miR-525-5p	
QM045_02	G18	HmiRQP0616	MIMAT0004780	hsa-miR-532-3p	



Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_02	G19	HmiRQP0615	MIMAT0002888	hsa-miR-532-5p	
QM045_02	G20	HmiRQP3008	MIMAT0022705	hsa-miR-539-3p	
QM045_02	G21	HmiRQP0621	MIMAT0003389	hsa-miR-542-3p	
QM045_02	G22	HmiRQP0622	MIMAT0004954	hsa-miR-543	
QM045_02	G23	HmiRQP0626	MIMAT0003251	hsa-miR-548a-3p	
QM045_02	G24	HmiRQP0627	MIMAT0004803	hsa-miR-548a-5p	
QM045_02	H01	HmiRQP2631	MIMAT0022266	hsa-miR-548ar-3p	
QM045_02	H02	HmiRQP0630	MIMAT0004806	hsa-miR-548c-5p	
QM045_02	H03	HmiRQP2015	MIMAT0018354	hsa-miR-548y	
QM045_02	H04	HmiRQP0665	MIMAT0003230	hsa-miR-566	
QM045_02	H05	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM045_02	H06	HmiRQP0674	MIMAT0004795	hsa-miR-574-5p	
QM045_02	H07	HmiRQP0675	MIMAT0003240	hsa-miR-575	
QM045_02	H08	HmiRQP0678	MIMAT0003242	hsa-miR-577	
QM045_02	H09	HmiRQP2753	MIMAT0023252	hsa-miR-5787	
QM045_02	H10	HmiRQP0680	MIMAT0003244	hsa-miR-579-3p	
QM045_02	H11	HmiRQP0684	MIMAT0003247	hsa-miR-582-5p	
QM045_02	H12	HmiRQP0686	MIMAT0003249	hsa-miR-584-5p	
QM045_02	H13	HmiRQP0703	MIMAT0003267	hsa-miR-599	
QM045_02	H14	HmiRQP2775	MIMAT0023714	hsa-miR-6089	
QM045_02	H15	HmiRQP0716	MIMAT0003280	hsa-miR-612	
QM045_02	H16	HmiRQP2779	MIMAT0024599	hsa-miR-6126	
QM045_02	H17	HmiRQP0720	MIMAT0003283	hsa-miR-615-3p	
QM045_02	H18	HmiRQP3265	MIMAT0026622	hsa-miR-619-5p	
QM045_02	H19	HmiRQP0734	MIMAT0003295	hsa-miR-626	
QM045_02	H20	HmiRQP0740	MIMAT0003299	hsa-miR-630	
QM045_02	H21	HmiRQP0748	MIMAT0003308	hsa-miR-638	
QM045_02	H22	HmiRQP3010	MIMAT0020924	hsa-miR-642a-3p	
QM045_02	H23	HmiRQP0760	MIMAT0003320	hsa-miR-650	
QM045_02	H24	HmiRQP0762	MIMAT0003322	hsa-miR-652-3p	
QM045_02	I01	HmiRQP3268	MIMAT0022709	hsa-miR-652-5p	
QM045_02	I02	HmiRQP0764	MIMAT0004814	hsa-miR-654-3p	
QM045_02	I03	HmiRQP0769	MIMAT0003336	hsa-miR-658	
QM045_02	I04	HmiRQP0778	MIMAT0004952	hsa-miR-665	
QM045_02	I05	HmiRQP0780	MIMAT0003880	hsa-miR-671-5p	
QM045_02	I06	HmiRQP3596	MIMAT0027400	hsa-miR-6750-5p	
QM045_02	I07	HmiRQP3599	MIMAT0027403	hsa-miR-6751-3p	
QM045_02	I08	HmiRQP3603	MIMAT0027407	hsa-miR-6753-3p	
QM045_02	I09	HmiRQP1310	MIMAT0006790	hsa-miR-675-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_02	I10	HmiRQP0782	MIMAT0004284	hsa-miR-675-5p	
QM045_02	I11	HmiRQP3622	MIMAT0027426	hsa-miR-6763-5p	
QM045_02	I12	HmiRQP2007	MIMAT0018204	hsa-miR-676-3p	
QM045_02	I13	HmiRQP3689	MIMAT0027493	hsa-miR-6796-3p	
QM045_02	I14	HmiRQP3703	MIMAT0027507	hsa-miR-6803-3p	
QM045_02	I15	HmiRQP0784	MIMAT0004927	hsa-miR-708-3p	
QM045_02	I16	HmiRQP0789	MIMAT0004946	hsa-miR-744-3p	
QM045_02	I17	HmiRQP0790	MIMAT0004945	hsa-miR-744-5p	
QM045_02	I18	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM045_02	I19	HmiRQP0792	MIMAT0004957	hsa-miR-760	
QM045_02	I20	HmiRQP1158	MIMAT0010364	hsa-miR-761	
QM045_02	I21	HmiRQP0796	MIMAT0003882	hsa-miR-767-5p	
QM045_02	I22	HmiRQP4536	MIMAT0030019	hsa-miR-7704	
QM045_02	I23	HmiRQP0801	MIMAT0003948	hsa-miR-770-5p	
QM045_02	I24	HmiRQP4538	MIMAT0030021	hsa-miR-7706	
QM045_02	J01	HmiRQP4563	MIMAT0031178	hsa-miR-7975	
QM045_02	J02	HmiRQP0802	MIMAT0004185	hsa-miR-802	
QM045_02	J03	HmiRQP3342	MIMAT0022717	hsa-miR-873-3p	
QM045_02	J04	HmiRQP0804	MIMAT0004911	hsa-miR-874-3p	
QM045_02	J05	HmiRQP3334	MIMAT0026718	hsa-miR-874-5p	
QM045_02	J06	HmiRQP0809	MIMAT0004949	hsa-miR-877-5p	
QM045_02	J07	HmiRQP0812	MIMAT0004947	hsa-miR-885-5p	
QM045_02	J08	HmiRQP0817	MIMAT0004916	hsa-miR-888-5p	
QM045_02	J09	HmiRQP0831	MIMAT0004507	hsa-miR-92a-1-5p	
QM045_02	J10	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM045_02	J11	HmiRQP0835	MIMAT0003218	hsa-miR-92b-3p	
QM045_02	J12	HmiRQP0834	MIMAT0004792	hsa-miR-92b-5p	
QM045_02	J13	HmiRQP0838	MIMAT0004976	hsa-miR-933	
QM045_02	J14	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM045_02	J15	HmiRQP0839	MIMAT0004977	hsa-miR-934	
QM045_02	J16	HmiRQP0840	MIMAT0004978	hsa-miR-935	
QM045_02	J17	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM045_02	J18	HmiRQP0841	MIMAT0004979	hsa-miR-936	
QM045_02	J19	HmiRQP0842	MIMAT0004980	hsa-miR-937-3p	
QM045_02	J20	HmiRQP0843	MIMAT0004981	hsa-miR-938	
QM045_02	J21	HmiRQP3346	MIMAT0022939	hsa-miR-939-3p	
QM045_02	J22	HmiRQP0844	MIMAT0004982	hsa-miR-939-5p	
QM045_02	J23	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM045_02	J24	HmiRQP0845	MIMAT0004983	hsa-miR-940	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM045_02	K01	HmiRQP0846	MIMAT0004984	hsa-miR-941	
QM045_02	K02	HmiRQP0849	MIMAT0004987	hsa-miR-944	
QM045_02	K03	HmiRQP3017	MIMAT0026473	hsa-miR-95-5p	
QM045_02	K04	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM045_02	K05	HmiRQP0851	MIMAT0004510	hsa-miR-96-3p	
QM045_02	K06	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM045_02	K07	HmiRQP0853	MIMAT0000096	hsa-miR-98-5p	
QM045_02	K08	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM045_02	K09	HmiRQP0856	MIMAT0004678	hsa-miR-99b-3p	
QM045_02	K10	HmiRQP0857	MIMAT0000689	hsa-miR-99b-5p	
QM045_02	L01	NC	NC	NC	
QM045_02	L02	NC	NC	NC	
QM045_02	L03	NC	NC	NC	
QM045_02	L04	NC	NC	NC	
QM045_02	L05	HmiRQP9001	NR_002752	HK1	
QM045_02	L06	HmiRQP9001	NR_002752	HK1	
QM045_02	L07	HmiRQP9011	NR_002750	HK2	
QM045_02	L08	HmiRQP9011	NR_002750	HK2	
QM045_02	L09	HmiRQP9021	NR_002745	HK3	
QM045_02	L10	HmiRQP9021	NR_002745	HK3	
QM045_02	L11	HmiRQP9051	NR_002746	HK4	
QM045_02	L12	HmiRQP9051	NR_002746	HK4	
QM045_02	L13	HmiRQP9061	NR_002744	HK5	
QM045_02	L14	HmiRQP9061	NR_002744	HK5	
QM045_02	L15	HmiRQP9071	NR_002450	HK6	
QM045_02	L16	HmiRQP9071	NR_002450	HK6	
QM045_02	L17	RT	RT	RT	
QM045_02	L18	RT	RT	RT	
QM045_02	L19	RT	RT	RT	
QM045_02	L20	RT	RT	RT	
QM045_02	L21	PCR	PCR	PCR	
QM045_02	L22	PCR	PCR	PCR	
QM045_02	L23	PCR	PCR	PCR	
QM045_02	L24	PCR	PCR	PCR	