

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM038-01	A01	MmiRQP0029	MIMAT0000386	mmu-miR-106b-5p	
QM038-01	A02	MmiRQP0098	MIMAT0000137	mmu-miR-126-5p	
QM038-01	A03	MmiRQP0125	MIMAT0000140	mmu-miR-128-3p	
QM038-01	A04	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM038-01	A05	MmiRQP0190	MIMAT0000156	mmu-miR-144-3p	
QM038-01	A06	MmiRQP0223	MIMAT0000526	mmu-miR-15a-5p	
QM038-01	A07	MmiRQP0225	MIMAT0000124	mmu-miR-15b-5p	
QM038-01	A08	MmiRQP0227	MIMAT0000527	mmu-miR-16-5p	
QM038-01	A09	MmiRQP0248	MIMAT0000215	mmu-miR-186-5p	
QM038-01	A10	MmiRQP0283	MIMAT0000225	mmu-miR-195-5p	
QM038-01	A11	MmiRQP0302	MIMAT0000657	mmu-miR-200c-3p	
QM038-01	A12	MmiRQP0338	MIMAT0000669	mmu-miR-221-3p	
QM038-01	B01	MmiRQP0339	MIMAT0000670	mmu-miR-222-3p	
QM038-01	B02	MmiRQP0354	MIMAT0000533	mmu-miR-26a-5p	
QM038-01	B03	MmiRQP0357	MIMAT0000534	mmu-miR-26b-5p	
QM038-01	B04	MmiRQP0373	MIMAT0000127	mmu-miR-29b-3p	
QM038-01	B05	MmiRQP0393	MIMAT0000130	mmu-miR-30b-5p	
QM038-01	B06	MmiRQP0396	MIMAT0000514	mmu-miR-30c-5p	
QM038-01	B07	MmiRQP0434	MIMAT0004651	mmu-miR-340-5p	
QM038-01	B08	MmiRQP0454	MIMAT0003186	mmu-miR-369-3p	
QM038-01	B09	MmiRQP0464	MIMAT0003727	mmu-miR-374-5p	
QM038-01	B10	MmiRQP0485	MIMAT0001091	mmu-miR-410-3p	
QM038-01	B11	MmiRQP0537	MIMAT0003456	mmu-miR-495-3p	
QM038-01	B12	MmiRQP0694	MIMAT0004896	mmu-miR-590-3p	
QM038-01	C01	MmiRQP0825	MIMAT0000142	mmu-miR-9-5p	
QM038-01	C02	MmiRQP0837	MIMAT0000540	mmu-miR-93-5p	
QM038-01	C03	MmiRQP0869	MIMAT0005850	mmu-miR-1192	
QM038-01	C04	MmiRQP0890	MIMAT0000165	mmu-miR-155-5p	
QM038-01	C05	MmiRQP0895	MIMAT0000211	mmu-miR-182-5p	
QM038-01	C06	MmiRQP0947	MIMAT0000372	mmu-miR-294-3p	
QM038-01	C07	MmiRQP0949	MIMAT0000373	mmu-miR-295-3p	
QM038-01	C08	MmiRQP0968	MIMAT0004640	mmu-miR-325-3p	
QM038-01	C09	MmiRQP0982	MIMAT0000605	mmu-miR-350-3p	
QM038-01	C10	MmiRQP0999	MIMAT0001537	mmu-miR-429-3p	
QM038-01	C11	MmiRQP1023	MIMAT0004930	mmu-miR-466d-5p	
QM038-01	C12	MmiRQP1026	MIMAT0004882	mmu-miR-466f-3p	
QM038-01	D01	MmiRQP1058	MIMAT0003453	mmu-miR-497-5p	
QM038-01	D02	MmiRQP1089	MIMAT0005839	mmu-miR-669f-3p	
QM038-01	D03	MmiRQP1123	MIMAT0003474	mmu-miR-694	
QM038-01	D04	MmiRQP1140	MIMAT0003502	mmu-miR-712-5p	
QM038-01	D05	MmiRQP1149	MIMAT0003515	mmu-miR-721	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM038-01	D06	MmiRQP1153	MIMAT0004238	mmu-miR-743a-3p	
QM038-01	D07	MmiRQP1154	MIMAT0004840	mmu-miR-743b-3p	
QM038-01	D08	MmiRQP1171	MIMAT0004855	mmu-miR-876-3p	
QM038-01	D09	MmiRQP1226	MIMAT0000548	mmu-miR-322-5p	
QM038-01	D10	MmiRQP1259	MIMAT0004745	mmu-miR-384-5p	
QM038-01	D11	MmiRQP0007	MIMAT0000383	mmu-let-7d-5p	
QM038-01	D12	MmiRQP0015	MIMAT0000121	mmu-let-7g-5p	
QM038-01	E01	MmiRQP0017	MIMAT0000122	mmu-let-7i-5p	
QM038-01	E02	MmiRQP0156	MIMAT0000141	mmu-miR-130a-3p	
QM038-01	E03	MmiRQP0159	MIMAT0000387	mmu-miR-130b-3p	
QM038-01	E04	MmiRQP0312	MIMAT0000529	mmu-miR-20a-5p	
QM038-01	E05	MmiRQP0314	MIMAT0003187	mmu-miR-20b-5p	
QM038-01	E06	MmiRQP0387	MIMAT0003377	mmu-miR-302d-3p	
QM038-01	E07	MmiRQP0853	MIMAT0000545	mmu-miR-98-5p	
QM038-01	E08	MmiRQP1092	MIMAT0005842	mmu-miR-669h-3p	
QM038-01	E09	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM038-01	E10	MmiRQP0004	MIMAT0000522	mmu-let-7b-5p	
QM038-01	E11	MmiRQP0006	MIMAT0000523	mmu-let-7c-5p	
QM038-01	E12	MmiRQP0010	MIMAT0000524	mmu-let-7e-5p	
QM038-01	F01	MmiRQP0012	MIMAT0000525	mmu-let-7f-5p	
QM038-01	F02	MmiRQP0170	MIMAT0000147	mmu-miR-135a-5p	
QM038-01	F03	MmiRQP0230	MIMAT0000649	mmu-miR-17-5p	
QM038-01	F04	MmiRQP0232	MIMAT0000210	mmu-miR-181a-5p	
QM038-01	F05	MmiRQP0234	MIMAT0000673	mmu-miR-181b-5p	
QM038-01	F06	MmiRQP0235	MIMAT0000674	mmu-miR-181c-5p	
QM038-01	F07	MmiRQP0237	MIMAT0004324	mmu-miR-181d-5p	
QM038-01	F08	MmiRQP0293	MIMAT0000651	mmu-miR-19a-3p	
QM038-01	F09	MmiRQP0295	MIMAT0000513	mmu-miR-19b-3p	
QM038-01	F10	MmiRQP0344	MIMAT0000532	mmu-miR-23a-3p	
QM038-01	F11	MmiRQP0346	MIMAT0000125	mmu-miR-23b-3p	
QM038-01	F12	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM038-01	G01	MmiRQP0361	MIMAT0000126	mmu-miR-27b-3p	
QM038-01	G02	MmiRQP0371	MIMAT0000535	mmu-miR-29a-3p	
QM038-01	G03	MmiRQP0375	MIMAT0000536	mmu-miR-29c-3p	
QM038-01	G04	MmiRQP0378	MIMAT0000379	mmu-miR-301a-3p	
QM038-01	G05	MmiRQP0391	MIMAT0000128	mmu-miR-30a-5p	
QM038-01	G06	MmiRQP0398	MIMAT0000515	mmu-miR-30d-5p	
QM038-01	G07	MmiRQP0400	MIMAT0000248	mmu-miR-30e-5p	
QM038-01	G08	MmiRQP0862	MIMAT0000385	mmu-miR-106a-5p	
QM038-01	G09	MmiRQP0941	MIMAT0000368	mmu-miR-291a-3p	
QM038-01	G10	MmiRQP0959	MIMAT0004186	mmu-miR-301b-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM038-01	G11	MmiRQP1019	MIMAT0005453	mmu-miR-466b-3-3p	
QM038-01	G12	MmiRQP1022	MIMAT0004931	mmu-miR-466d-3p	
QM038-01	H01	NC	NC	NC	
QM038-01	H02	NC	NC	NC	
QM038-01	H03	MmiRQP9002	NR_003027	MK1	
QM038-01	H04	MmiRQP9012	NR_028128	MK2	
QM038-01	H05	MmiRQP9022	NR_028554	MK3	
QM038-01	H06	MmiRQP9042	NR_028550	MK4	
QM038-01	H07	MmiRQP9052	NR_028541	MK5	
QM038-01	H08	MmiRQP9062	NR_028091	MK6	
QM038-01	H09	RT	RT	RT	
QM038-01	H10	RT	RT	RT	
QM038-01	H11	PCR	PCR	PCR	
QM038-01	H12	PCR	PCR	PCR	