

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM024-01	A01	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM024-01	A02	MmiRQP0003	MIMAT0004621	mmu-let-7b-3p	
QM024-01	A03	MmiRQP0004	MIMAT0000522	mmu-let-7b-5p	
QM024-01	A04	MmiRQP0006	MIMAT0000523	mmu-let-7c-5p	
QM024-01	A05	MmiRQP0008	MIMAT0000384	mmu-let-7d-3p	
QM024-01	A06	MmiRQP0007	MIMAT0000383	mmu-let-7d-5p	
QM024-01	A07	MmiRQP0010	MIMAT0000524	mmu-let-7e-5p	
QM024-01	A08	MmiRQP0012	MIMAT0000525	mmu-let-7f-5p	
QM024-01	A09	MmiRQP0015	MIMAT0000121	mmu-let-7g-5p	
QM024-01	A10	MmiRQP0016	MIMAT0004520	mmu-let-7i-3p	
QM024-01	A11	MmiRQP0017	MIMAT0000122	mmu-let-7i-5p	
QM024-01	A12	MmiRQP0018	MIMAT0000655	mmu-miR-100-5p	
QM024-01	B01	MmiRQP0056	MIMAT0000246	mmu-miR-122-5p	
QM024-01	B02	MmiRQP0074	MIMAT0000134	mmu-miR-124-3p	
QM024-01	B03	MmiRQP0096	MIMAT0000136	mmu-miR-125b-5p	
QM024-01	B04	MmiRQP0111	MIMAT0000139	mmu-miR-127-3p	
QM024-01	B05	MmiRQP0110	MIMAT0004530	mmu-miR-127-5p	
QM024-01	B06	MmiRQP0125	MIMAT0000140	mmu-miR-128-3p	
QM024-01	B07	MmiRQP0139	MIMAT0000544	mmu-miR-129-2-3p	
QM024-01	B08	MmiRQP0137	MIMAT0000209	mmu-miR-129-5p	
QM024-01	B09	MmiRQP0170	MIMAT0000147	mmu-miR-135a-5p	
QM024-01	B10	MmiRQP0172	MIMAT0000612	mmu-miR-135b-5p	
QM024-01	B11	MmiRQP0175	MIMAT0000149	mmu-miR-137-3p	
QM024-01	B12	MmiRQP0180	MIMAT0000656	mmu-miR-139-5p	
QM024-01	C01	MmiRQP0182	MIMAT0000152	mmu-miR-140-3p	
QM024-01	C02	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM024-01	C03	MmiRQP0184	MIMAT0000153	mmu-miR-141-3p	
QM024-01	C04	MmiRQP0885	MIMAT0004533	mmu-miR-141-5p	
QM024-01	C05	MmiRQP0188	MIMAT0000247	mmu-miR-143-3p	
QM024-01	C06	MmiRQP0190	MIMAT0000156	mmu-miR-144-3p	
QM024-01	C07	MmiRQP0208	MIMAT0000159	mmu-miR-149-5p	
QM024-01	C08	MmiRQP0219	MIMAT0000164	mmu-miR-154-5p	
QM024-01	C09	MmiRQP0223	MIMAT0000526	mmu-miR-15a-5p	
QM024-01	C10	MmiRQP0225	MIMAT0000124	mmu-miR-15b-5p	
QM024-01	C11	MmiRQP0232	MIMAT0000210	mmu-miR-181a-5p	
QM024-01	C12	MmiRQP0243	MIMAT0004539	mmu-miR-183-3p	
QM024-01	D01	MmiRQP0244	MIMAT0000212	mmu-miR-183-5p	
QM024-01	D02	MmiRQP0245	MIMAT0000213	mmu-miR-184-3p	
QM024-01	D03	MmiRQP0247	MIMAT0000214	mmu-miR-185-5p	
QM024-01	D04	MmiRQP0258	MIMAT0000220	mmu-miR-190-5p	
QM024-01	D05	MmiRQP0263	MIMAT0000221	mmu-miR-191-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM024-01	D06	MmiRQP0276	MIMAT0000223	mmu-miR-193-3p	
QM024-01	D07	MmiRQP0284	MIMAT0000518	mmu-miR-196a-5p	
QM024-01	D08	MmiRQP0290	MIMAT0000229	mmu-miR-199a-5p	
QM024-01	D09	MmiRQP0293	MIMAT0000651	mmu-miR-19a-3p	
QM024-01	D10	MmiRQP0298	MIMAT0000519	mmu-miR-200a-3p	
QM024-01	D11	MmiRQP0300	MIMAT0000233	mmu-miR-200b-3p	
QM024-01	D12	MmiRQP0299	MIMAT0004545	mmu-miR-200b-5p	
QM024-01	E01	MmiRQP0302	MIMAT0000657	mmu-miR-200c-3p	
QM024-01	E02	MmiRQP0301	MIMAT0004663	mmu-miR-200c-5p	
QM024-01	E03	MmiRQP0306	MIMAT0000237	mmu-miR-204-5p	
QM024-01	E04	MmiRQP0317	MIMAT0000658	mmu-miR-210-3p	
QM024-01	E05	MmiRQP0929	MIMAT0004628	mmu-miR-21-3p	
QM024-01	E06	MmiRQP0316	MIMAT0000530	mmu-miR-21-5p	
QM024-01	E07	MmiRQP0324	MIMAT0003729	mmu-miR-216b-5p	
QM024-01	E08	MmiRQP0330	MIMAT0000664	mmu-miR-219-5p	
QM024-01	E09	MmiRQP0338	MIMAT0000669	mmu-miR-221-3p	
QM024-01	E10	MmiRQP0339	MIMAT0000670	mmu-miR-222-3p	
QM024-01	E11	MmiRQP0346	MIMAT0000125	mmu-miR-23b-3p	
QM024-01	E12	MmiRQP0937	MIMAT0000218	mmu-miR-24-1-5p	
QM024-01	F01	MmiRQP0349	MIMAT0000219	mmu-miR-24-3p	
QM024-01	F02	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM024-01	F03	MmiRQP0358	MIMAT0004633	mmu-miR-27a-5p	
QM024-01	F04	MmiRQP0362	MIMAT0000653	mmu-miR-28-5p	
QM024-01	F05	MmiRQP0373	MIMAT0000127	mmu-miR-29b-3p	
QM024-01	F06	MmiRQP0381	MIMAT0000380	mmu-miR-302a-3p	
QM024-01	F07	MmiRQP0387	MIMAT0003377	mmu-miR-302d-3p	
QM024-01	F08	MmiRQP0393	MIMAT0000130	mmu-miR-30b-5p	
QM024-01	F09	MmiRQP0965	MIMAT0004634	mmu-miR-31-3p	
QM024-01	F10	MmiRQP1224	MIMAT0000538	mmu-miR-31-5p	
QM024-01	F11	MmiRQP0410	MIMAT0000551	mmu-miR-323-3p	
QM024-01	F12	MmiRQP0409	MIMAT0004638	mmu-miR-323-5p	
QM024-01	G01	MmiRQP0412	MIMAT0000555	mmu-miR-324-5p	
QM024-01	G02	MmiRQP0420	MIMAT0000571	mmu-miR-331-3p	
QM024-01	G03	MmiRQP0419	MIMAT0004643	mmu-miR-331-5p	
QM024-01	G04	MmiRQP0426	MIMAT0000582	mmu-miR-338-3p	
QM024-01	G05	MmiRQP0440	MIMAT0000542	mmu-miR-34a-5p	
QM024-01	G06	MmiRQP0444	MIMAT0000381	mmu-miR-34c-5p	
QM024-01	G07	MmiRQP0464	MIMAT0003727	mmu-miR-374-5p	
QM024-01	G08	MmiRQP0615	MIMAT0002889	mmu-miR-532-5p	
QM024-01	G09	MmiRQP0762	MIMAT0003711	mmu-miR-652-3p	
QM024-01	G10	MmiRQP0825	MIMAT0000142	mmu-miR-9-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM024-01	G11	MmiRQP0852	MIMAT0000541	mmu-miR-96-5p	
QM024-01	G12	MmiRQP0853	MIMAT0000545	mmu-miR-98-5p	
QM024-01	H01	NC	NC	NC	
QM024-01	H02	NC	NC	NC	
QM024-01	H03	MmiRQP9002	NR_003027	MK1	
QM024-01	H04	MmiRQP9012	NR_028128	MK2	
QM024-01	H05	MmiRQP9022	NR_028554	MK3	
QM024-01	H06	MmiRQP9042	NR_028550	MK4	
QM024-01	H07	MmiRQP9052	NR_028541	MK5	
QM024-01	H08	MmiRQP9062	NR_028091	MK6	
QM024-01	H09	RT	RT	RT	
QM024-01	H10	RT	RT	RT	
QM024-01	H11	PCR	PCR	PCR	
QM024-01	H12	PCR	PCR	PCR	