

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM026-01	A01	MmiRQP0029	MIMAT0000386	mmu-miR-106b-5p	
QM026-01	A02	MmiRQP0096	MIMAT0000136	mmu-miR-125b-5p	
QM026-01	A03	MmiRQP0099	MIMAT0000138	mmu-miR-126-3p	
QM026-01	A04	MmiRQP0168	MIMAT0000146	mmu-miR-134-5p	
QM026-01	A05	MmiRQP0175	MIMAT0000149	mmu-miR-137-3p	
QM026-01	A06	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM026-01	A07	MmiRQP0185	MIMAT0000154	mmu-miR-142-5p	
QM026-01	A08	MmiRQP0188	MIMAT0000247	mmu-miR-143-3p	
QM026-01	A09	MmiRQP0192	MIMAT0000157	mmu-miR-145-5p	
QM026-01	A10	MmiRQP0219	MIMAT0000164	mmu-miR-154-5p	
QM026-01	A11	MmiRQP0223	MIMAT0000526	mmu-miR-15a-5p	
QM026-01	A12	MmiRQP0227	MIMAT0000527	mmu-miR-16-5p	
QM026-01	B01	MmiRQP0245	MIMAT0000213	mmu-miR-184-3p	
QM026-01	B02	MmiRQP0258	MIMAT0000220	mmu-miR-190-5p	
QM026-01	B03	MmiRQP0274	MIMAT0000517	mmu-miR-192-5p	
QM026-01	B04	MmiRQP0283	MIMAT0000225	mmu-miR-195-5p	
QM026-01	B05	MmiRQP0290	MIMAT0000229	mmu-miR-199a-5p	
QM026-01	B06	MmiRQP0300	MIMAT0000233	mmu-miR-200b-3p	
QM026-01	B07	MmiRQP0302	MIMAT0000657	mmu-miR-200c-3p	
QM026-01	B08	MmiRQP0305	MIMAT0000236	mmu-miR-203-3p	
QM026-01	B09	MmiRQP0307	MIMAT0000238	mmu-miR-205-5p	
QM026-01	B10	MmiRQP0308	MIMAT0000239	mmu-miR-206-3p	
QM026-01	B11	MmiRQP0316	MIMAT0000530	mmu-miR-21-5p	
QM026-01	B12	MmiRQP0320	MIMAT0000661	mmu-miR-214-3p	
QM026-01	C01	MmiRQP0332	MIMAT0000531	mmu-miR-22-3p	
QM026-01	C02	MmiRQP0338	MIMAT0000669	mmu-miR-221-3p	
QM026-01	C03	MmiRQP0339	MIMAT0000670	mmu-miR-222-3p	
QM026-01	C04	MmiRQP0342	MIMAT0000665	mmu-miR-223-3p	
QM026-01	C05	MmiRQP0354	MIMAT0000533	mmu-miR-26a-5p	
QM026-01	C06	MmiRQP0421	MIMAT0000766	mmu-miR-335-5p	
QM026-01	C07	MmiRQP0440	MIMAT0000542	mmu-miR-34a-5p	
QM026-01	C08	MmiRQP0451	MIMAT0000711	mmu-miR-365-3p	
QM026-01	C09	MmiRQP0466	MIMAT0000739	mmu-miR-375-3p	
QM026-01	C10	MmiRQP0472	MIMAT0000741	mmu-miR-377-3p	
QM026-01	C11	MmiRQP0476	MIMAT0000743	mmu-miR-379-5p	
QM026-01	C12	MmiRQP0485	MIMAT0001091	mmu-miR-410-3p	
QM026-01	D01	MmiRQP0503	MIMAT0001533	mmu-miR-448-3p	
QM026-01	D02	MmiRQP0525	MIMAT0003184	mmu-miR-487b-3p	
QM026-01	D03	MmiRQP0537	MIMAT0003456	mmu-miR-495-3p	
QM026-01	D04	MmiRQP0825	MIMAT0000142	mmu-miR-9-5p	
QM026-01	D05	MmiRQP0837	MIMAT0000540	mmu-miR-93-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM026-01	D06	MmiRQP0861	MIMAT0004856	mmu-miR-105	
QM026-01	D07	MmiRQP0890	MIMAT0000165	mmu-miR-155-5p	
QM026-01	D08	MmiRQP0895	MIMAT0000211	mmu-miR-182-5p	
QM026-01	D09	MmiRQP0936	MIMAT0000671	mmu-miR-224-5p	
QM026-01	D10	MmiRQP0968	MIMAT0004640	mmu-miR-325-3p	
QM026-01	D11	MmiRQP0978	MIMAT0000597	mmu-miR-346-5p	
QM026-01	D12	MmiRQP0987	MIMAT0000740	mmu-miR-376a-3p	
QM026-01	E01	MmiRQP0999	MIMAT0001537	mmu-miR-429-3p	
QM026-01	E02	MmiRQP1055	MIMAT0003450	mmu-miR-488-3p	
QM026-01	E03	MmiRQP0007	MIMAT0000383	mmu-let-7d-5p	
QM026-01	E04	MmiRQP0017	MIMAT0000122	mmu-let-7i-5p	
QM026-01	E05	MmiRQP0184	MIMAT0000153	mmu-miR-141-3p	
QM026-01	E06	MmiRQP0196	MIMAT0000158	mmu-miR-146a-5p	
QM026-01	E07	MmiRQP0298	MIMAT0000519	mmu-miR-200a-3p	
QM026-01	E08	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM026-01	E09	MmiRQP0004	MIMAT0000522	mmu-let-7b-5p	
QM026-01	E10	MmiRQP0006	MIMAT0000523	mmu-let-7c-5p	
QM026-01	E11	MmiRQP0018	MIMAT0000655	mmu-miR-100-5p	
QM026-01	E12	MmiRQP0021	MIMAT0000133	mmu-miR-101a-3p	
QM026-01	F01	MmiRQP0034	MIMAT0000208	mmu-miR-10b-5p	
QM026-01	F02	MmiRQP0166	MIMAT0000145	mmu-miR-133a-3p	
QM026-01	F03	MmiRQP0306	MIMAT0000237	mmu-miR-204-5p	
QM026-01	F04	MmiRQP0344	MIMAT0000532	mmu-miR-23a-3p	
QM026-01	F05	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM026-01	F06	MmiRQP0371	MIMAT0000535	mmu-miR-29a-3p	
QM026-01	F07	MmiRQP0375	MIMAT0000536	mmu-miR-29c-3p	
QM026-01	F08	MmiRQP0391	MIMAT0000128	mmu-miR-30a-5p	
QM026-01	F09	MmiRQP0400	MIMAT0000248	mmu-miR-30e-5p	
QM026-01	F10	MmiRQP0854	MIMAT0000131	mmu-miR-99a-5p	
QM026-01	F11	MmiRQP0930	MIMAT0000668	mmu-miR-211-5p	
QM026-01	F12	MmiRQP0289	MIMAT0004667	mmu-miR-199b-3p	mmu-miR-199a-3p
QM026-01	G01	MmiRQP0218	MIMAT0004537	mmu-miR-154-3p	
QM026-01	G02	MmiRQP2568	MIMAT0017022	mmu-miR-34a-3p	
QM026-01	G03	MmiRQP0929	MIMAT0004628	mmu-miR-21-3p	
QM026-01	G04	MmiRQP0028	MIMAT0004582	mmu-miR-106b-3p	
QM026-01	G05	MmiRQP0182	MIMAT0000152	mmu-miR-140-3p	
QM026-01	G06	MmiRQP0390	MIMAT0000129	mmu-miR-30a-3p	
QM026-01	G07	MmiRQP0863	MIMAT0004538	mmu-miR-10b-3p	
QM026-01	G08	MmiRQP0883	MIMAT0004668	mmu-miR-138-1-3p	
QM026-01	G09	MmiRQP0925	MIMAT0004547	mmu-miR-203-5p	
QM026-01	G10	MmiRQP0961	MIMAT0003373	mmu-miR-302b-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM026-01	G11	MmiRQP0963	MIMAT0003375	mmu-miR-302c-5p	
QM026-01	G12	MmiRQP1228	MIMAT0000558	mmu-miR-325-5p	
QM026-01	H01	NC	NC	NC	
QM026-01	H02	NC	NC	NC	
QM026-01	H03	MmiRQP9002	NR_003027	MK1	
QM026-01	H04	MmiRQP9012	NR_028128	MK2	
QM026-01	H05	MmiRQP9022	NR_028554	MK3	
QM026-01	H06	MmiRQP9042	NR_028550	MK4	
QM026-01	H07	MmiRQP9052	NR_028541	MK5	
QM026-01	H08	MmiRQP9062	NR_028091	MK6	
QM026-01	H09	RT	RT	RT	
QM026-01	H10	RT	RT	RT	
QM026-01	H11	PCR	PCR	PCR	
QM026-01	H12	PCR	PCR	PCR	