

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM040-01	A01	MmiRQP0096	MIMAT0000136	mmu-miR-125b-5p	
QM040-01	A02	MmiRQP0099	MIMAT0000138	mmu-miR-126-3p	
QM040-01	A03	MmiRQP0137	MIMAT0000209	mmu-miR-129-5p	
QM040-01	A04	MmiRQP0161	MIMAT0000144	mmu-miR-132-3p	
QM040-01	A05	MmiRQP0168	MIMAT0000146	mmu-miR-134-5p	
QM040-01	A06	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM040-01	A07	MmiRQP0185	MIMAT0000154	mmu-miR-142-5p	
QM040-01	A08	MmiRQP0186	MIMAT0000155	mmu-miR-142-3p	
QM040-01	A09	MmiRQP0192	MIMAT0000157	mmu-miR-145-5p	
QM040-01	A10	MmiRQP0208	MIMAT0000159	mmu-miR-149-5p	
QM040-01	A11	MmiRQP0210	MIMAT0000160	mmu-miR-150-5p	
QM040-01	A12	MmiRQP0223	MIMAT0000526	mmu-miR-15a-5p	
QM040-01	B01	MmiRQP0225	MIMAT0000124	mmu-miR-15b-5p	
QM040-01	B02	MmiRQP0227	MIMAT0000527	mmu-miR-16-5p	
QM040-01	B03	MmiRQP0244	MIMAT0000212	mmu-miR-183-5p	
QM040-01	B04	MmiRQP0245	MIMAT0000213	mmu-miR-184-3p	
QM040-01	B05	MmiRQP0247	MIMAT0000214	mmu-miR-185-5p	
QM040-01	B06	MmiRQP0248	MIMAT0000215	mmu-miR-186-5p	
QM040-01	B07	MmiRQP0263	MIMAT0000221	mmu-miR-191-5p	
QM040-01	B08	MmiRQP0280	MIMAT0000224	mmu-miR-194-5p	
QM040-01	B09	MmiRQP0283	MIMAT0000225	mmu-miR-195-5p	
QM040-01	B10	MmiRQP0305	MIMAT0000236	mmu-miR-203-3p	
QM040-01	B11	MmiRQP0307	MIMAT0000238	mmu-miR-205-5p	
QM040-01	B12	MmiRQP0316	MIMAT0000530	mmu-miR-21-5p	
QM040-01	C01	MmiRQP0317	MIMAT0000658	mmu-miR-210-3p	
QM040-01	C02	MmiRQP0320	MIMAT0000661	mmu-miR-214-3p	
QM040-01	C03	MmiRQP0342	MIMAT0000665	mmu-miR-223-3p	
QM040-01	C04	MmiRQP0354	MIMAT0000533	mmu-miR-26a-5p	
QM040-01	C05	MmiRQP0357	MIMAT0000534	mmu-miR-26b-5p	
QM040-01	C06	MmiRQP0362	MIMAT0000653	mmu-miR-28-5p	
QM040-01	C07	MmiRQP0373	MIMAT0000127	mmu-miR-29b-3p	
QM040-01	C08	MmiRQP0393	MIMAT0000130	mmu-miR-30b-5p	
QM040-01	C09	MmiRQP0396	MIMAT0000514	mmu-miR-30c-5p	
QM040-01	C10	MmiRQP0421	MIMAT0000766	mmu-miR-335-5p	
QM040-01	C11	MmiRQP0440	MIMAT0000542	mmu-miR-34a-5p	
QM040-01	C12	MmiRQP0449	MIMAT0000708	mmu-miR-363-3p	
QM040-01	D01	MmiRQP0456	MIMAT0001095	mmu-miR-370-3p	
QM040-01	D02	MmiRQP0484	MIMAT0001090	mmu-miR-409-3p	
QM040-01	D03	MmiRQP0509	MIMAT0001632	mmu-miR-451	
QM040-01	D04	MmiRQP0673	MIMAT0004894	mmu-miR-574-3p	
QM040-01	D05	MmiRQP0825	MIMAT0000142	mmu-miR-9-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM040-01	D06	MmiRQP0857	MIMAT0000132	mmu-miR-99b-5p	
QM040-01	D07	MmiRQP0861	MIMAT0004856	mmu-miR-105	
QM040-01	D08	MmiRQP0890	MIMAT0000165	mmu-miR-155-5p	
QM040-01	D09	MmiRQP0895	MIMAT0000211	mmu-miR-182-5p	
QM040-01	D10	MmiRQP0955	MIMAT0000376	mmu-miR-298-5p	
QM040-01	D11	MmiRQP0956	MIMAT0004577	mmu-miR-299-3p	
QM040-01	D12	MmiRQP0968	MIMAT0004640	mmu-miR-325-3p	
QM040-01	E01	MmiRQP0995	MIMAT0000748	mmu-miR-383-5p	
QM040-01	E02	MmiRQP1224	MIMAT0000538	mmu-miR-31-5p	
QM040-01	E03	MmiRQP0007	MIMAT0000383	mmu-let-7d-5p	
QM040-01	E04	MmiRQP0015	MIMAT0000121	mmu-let-7g-5p	
QM040-01	E05	MmiRQP0156	MIMAT0000141	mmu-miR-130a-3p	
QM040-01	E06	MmiRQP0196	MIMAT0000158	mmu-miR-146a-5p	
QM040-01	E07	MmiRQP0197	MIMAT0003475	mmu-miR-146b-5p	
QM040-01	E08	MmiRQP0204	MIMAT0000516	mmu-miR-148a-3p	
QM040-01	E09	MmiRQP0255	MIMAT0000528	mmu-miR-18a-5p	
QM040-01	E10	MmiRQP0298	MIMAT0000519	mmu-miR-200a-3p	
QM040-01	E11	MmiRQP0312	MIMAT0000529	mmu-miR-20a-5p	
QM040-01	E12	MmiRQP0314	MIMAT0003187	mmu-miR-20b-5p	
QM040-01	F01	MmiRQP0444	MIMAT0000381	mmu-miR-34c-5p	
QM040-01	F02	MmiRQP0853	MIMAT0000545	mmu-miR-98-5p	
QM040-01	F03	MmiRQP0908	MIMAT0004858	mmu-miR-18b-5p	
QM040-01	F04	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM040-01	F05	MmiRQP0006	MIMAT0000523	mmu-let-7c-5p	
QM040-01	F06	MmiRQP0010	MIMAT0000524	mmu-let-7e-5p	
QM040-01	F07	MmiRQP0012	MIMAT0000525	mmu-let-7f-5p	
QM040-01	F08	MmiRQP0170	MIMAT0000147	mmu-miR-135a-5p	
QM040-01	F09	MmiRQP0172	MIMAT0000612	mmu-miR-135b-5p	
QM040-01	F10	MmiRQP0232	MIMAT0000210	mmu-miR-181a-5p	
QM040-01	F11	MmiRQP0284	MIMAT0000518	mmu-miR-196a-5p	
QM040-01	F12	MmiRQP0293	MIMAT0000651	mmu-miR-19a-3p	
QM040-01	G01	MmiRQP0295	MIMAT0000513	mmu-miR-19b-3p	
QM040-01	G02	MmiRQP0346	MIMAT0000125	mmu-miR-23b-3p	
QM040-01	G03	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM040-01	G04	MmiRQP0361	MIMAT0000126	mmu-miR-27b-3p	
QM040-01	G05	MmiRQP0375	MIMAT0000536	mmu-miR-29c-3p	
QM040-01	G06	MmiRQP0400	MIMAT0000248	mmu-miR-30e-5p	
QM040-01	G07	MmiRQP0862	MIMAT0000385	mmu-miR-106a-5p	
QM040-01	G08	MmiRQP0369	MIMAT0000377	mmu-miR-299-5p	
QM040-01	G09	MmiRQP0929	MIMAT0004628	mmu-miR-21-3p	
QM040-01	G10	MmiRQP0824	MIMAT0000143	mmu-miR-9-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM040-01	G11	MmiRQP0883	MIMAT0004668	mmu-miR-138-1-3p	
QM040-01	G12	MmiRQP0965	MIMAT0004634	mmu-miR-31-3p	
QM040-01	H01	NC	NC	NC	
QM040-01	H02	NC	NC	NC	
QM040-01	H03	MmiRQP9002	NR_003027	MK1	
QM040-01	H04	MmiRQP9012	NR_028128	MK2	
QM040-01	H05	MmiRQP9022	NR_028554	MK3	
QM040-01	H06	MmiRQP9042	NR_028550	MK4	
QM040-01	H07	MmiRQP9052	NR_028541	MK5	
QM040-01	H08	MmiRQP9062	NR_028091	MK6	
QM040-01	H09	RT	RT	RT	
QM040-01	H10	RT	RT	RT	
QM040-01	H11	PCR	PCR	PCR	
QM040-01	H12	PCR	PCR	PCR	