

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM027-01	A01	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM027-01	A02	MmiRQP0006	MIMAT0000523	mmu-let-7c-5p	
QM027-01	A03	MmiRQP0007	MIMAT0000383	mmu-let-7d-5p	
QM027-01	A04	MmiRQP0018	MIMAT0000655	mmu-miR-100-5p	
QM027-01	A05	MmiRQP0021	MIMAT0000133	mmu-miR-101a-3p	
QM027-01	A06	MmiRQP0029	MIMAT0000386	mmu-miR-106b-5p	
QM027-01	A07	MmiRQP0032	MIMAT0000648	mmu-miR-10a-5p	
QM027-01	A08	MmiRQP0056	MIMAT0000246	mmu-miR-122-5p	
QM027-01	A09	MmiRQP0093	MIMAT0004528	mmu-miR-125a-3p	
QM027-01	A10	MmiRQP0096	MIMAT0000136	mmu-miR-125b-5p	
QM027-01	A11	MmiRQP0111	MIMAT0000139	mmu-miR-127-3p	
QM027-01	A12	MmiRQP0125	MIMAT0000140	mmu-miR-128-3p	
QM027-01	B01	MmiRQP0159	MIMAT0000387	mmu-miR-130b-3p	
QM027-01	B02	MmiRQP0161	MIMAT0000144	mmu-miR-132-3p	
QM027-01	B03	MmiRQP0166	MIMAT0000145	mmu-miR-133a-3p	
QM027-01	B04	MmiRQP0168	MIMAT0000146	mmu-miR-134-5p	
QM027-01	B05	MmiRQP0172	MIMAT0000612	mmu-miR-135b-5p	
QM027-01	B06	MmiRQP0175	MIMAT0000149	mmu-miR-137-3p	
QM027-01	B07	MmiRQP0180	MIMAT0000656	mmu-miR-139-5p	
QM027-01	B08	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM027-01	B09	MmiRQP0184	MIMAT0000153	mmu-miR-141-3p	
QM027-01	B10	MmiRQP0186	MIMAT0000155	mmu-miR-142-3p	
QM027-01	B11	MmiRQP0188	MIMAT0000247	mmu-miR-143-3p	
QM027-01	B12	MmiRQP0192	MIMAT0000157	mmu-miR-145-5p	
QM027-01	C01	MmiRQP0196	MIMAT0000158	mmu-miR-146a-5p	
QM027-01	C02	MmiRQP0204	MIMAT0000516	mmu-miR-148a-3p	
QM027-01	C03	MmiRQP0206	MIMAT0000580	mmu-miR-148b-3p	
QM027-01	C04	MmiRQP0208	MIMAT0000159	mmu-miR-149-5p	
QM027-01	C05	MmiRQP0210	MIMAT0000160	mmu-miR-150-5p	
QM027-01	C06	MmiRQP0214	MIMAT0000163	mmu-miR-153-3p	
QM027-01	C07	MmiRQP0219	MIMAT0000164	mmu-miR-154-5p	
QM027-01	C08	MmiRQP0223	MIMAT0000526	mmu-miR-15a-5p	
QM027-01	C09	MmiRQP0227	MIMAT0000527	mmu-miR-16-5p	
QM027-01	C10	MmiRQP0230	MIMAT0000649	mmu-miR-17-5p	
QM027-01	C11	MmiRQP0232	MIMAT0000210	mmu-miR-181a-5p	
QM027-01	C12	MmiRQP0244	MIMAT0000212	mmu-miR-183-5p	
QM027-01	D01	MmiRQP0247	MIMAT0000214	mmu-miR-185-5p	
QM027-01	D02	MmiRQP0248	MIMAT0000215	mmu-miR-186-5p	
QM027-01	D03	MmiRQP0255	MIMAT0000528	mmu-miR-18a-5p	
QM027-01	D04	MmiRQP0258	MIMAT0000220	mmu-miR-190-5p	
QM027-01	D05	MmiRQP0263	MIMAT0000221	mmu-miR-191-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM027-01	D06	MmiRQP0274	MIMAT0000517	mmu-miR-192-5p	
QM027-01	D07	MmiRQP0280	MIMAT0000224	mmu-miR-194-5p	
QM027-01	D08	MmiRQP0283	MIMAT0000225	mmu-miR-195-5p	
QM027-01	D09	MmiRQP0284	MIMAT0000518	mmu-miR-196a-5p	
QM027-01	D10	MmiRQP0290	MIMAT0000229	mmu-miR-199a-5p	
QM027-01	D11	MmiRQP0293	MIMAT0000651	mmu-miR-19a-3p	
QM027-01	D12	MmiRQP0295	MIMAT0000513	mmu-miR-19b-3p	
QM027-01	E01	MmiRQP0044	MIMAT0000123	mmu-miR-1a-3p	
QM027-01	E02	MmiRQP0298	MIMAT0000519	mmu-miR-200a-3p	
QM027-01	E03	MmiRQP0305	MIMAT0000236	mmu-miR-203-3p	
QM027-01	E04	MmiRQP0306	MIMAT0000237	mmu-miR-204-5p	
QM027-01	E05	MmiRQP0307	MIMAT0000238	mmu-miR-205-5p	
QM027-01	E06	MmiRQP0308	MIMAT0000239	mmu-miR-206-3p	
QM027-01	E07	MmiRQP0312	MIMAT0000529	mmu-miR-20a-5p	
QM027-01	E08	MmiRQP0317	MIMAT0000658	mmu-miR-210-3p	
QM027-01	E09	MmiRQP0320	MIMAT0000661	mmu-miR-214-3p	
QM027-01	E10	MmiRQP0316	MIMAT0000530	mmu-miR-21-5p	
QM027-01	E11	MmiRQP0324	MIMAT0003729	mmu-miR-216b-5p	
QM027-01	E12	MmiRQP0327	MIMAT0000663	mmu-miR-218-5p	
QM027-01	F01	MmiRQP0330	MIMAT0000664	mmu-miR-219-5p	
QM027-01	F02	MmiRQP0338	MIMAT0000669	mmu-miR-221-3p	
QM027-01	F03	MmiRQP0339	MIMAT0000670	mmu-miR-222-3p	
QM027-01	F04	MmiRQP0342	MIMAT0000665	mmu-miR-223-3p	
QM027-01	F05	MmiRQP0332	MIMAT0000531	mmu-miR-22-3p	
QM027-01	F06	MmiRQP0344	MIMAT0000532	mmu-miR-23a-3p	
QM027-01	F07	MmiRQP0346	MIMAT0000125	mmu-miR-23b-3p	
QM027-01	F08	MmiRQP0349	MIMAT0000219	mmu-miR-24-3p	
QM027-01	F09	MmiRQP0352	MIMAT0000652	mmu-miR-25-3p	
QM027-01	F10	MmiRQP0354	MIMAT0000533	mmu-miR-26a-5p	
QM027-01	F11	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM027-01	F12	MmiRQP0364	MIMAT0000374	mmu-miR-296-5p	
QM027-01	G01	MmiRQP0371	MIMAT0000535	mmu-miR-29a-3p	
QM027-01	G02	MmiRQP0378	MIMAT0000379	mmu-miR-301a-3p	
QM027-01	G03	MmiRQP0391	MIMAT0000128	mmu-miR-30a-5p	
QM027-01	G04	MmiRQP1224	MIMAT0000538	mmu-miR-31-5p	
QM027-01	G05	MmiRQP0404	MIMAT0000654	mmu-miR-32-5p	
QM027-01	G06	MmiRQP0440	MIMAT0000542	mmu-miR-34a-5p	
QM027-01	G07	MmiRQP0466	MIMAT0000739	mmu-miR-375-3p	
QM027-01	G08	MmiRQP0786	MIMAT0000677	mmu-miR-7a-5p	
QM027-01	G09	MmiRQP0837	MIMAT0000540	mmu-miR-93-5p	
QM027-01	G10	MmiRQP0825	MIMAT0000142	mmu-miR-9-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM027-01	G11	MmiRQP0852	MIMAT0000541	mmu-miR-96-5p	
QM027-01	G12	MmiRQP0854	MIMAT0000131	mmu-miR-99a-5p	
QM027-01	H01	NC	NC	NC	
QM027-01	H02	NC	NC	NC	
QM027-01	H03	MmiRQP9002	NR_003027	MK1	
QM027-01	H04	MmiRQP9012	NR_028128	MK2	
QM027-01	H05	MmiRQP9022	NR_028554	MK3	
QM027-01	H06	MmiRQP9042	NR_028550	MK4	
QM027-01	H07	MmiRQP9052	NR_028541	MK5	
QM027-01	H08	MmiRQP9062	NR_028091	MK6	
QM027-01	H09	RT	RT	RT	
QM027-01	H10	RT	RT	RT	
QM027-01	H11	PCR	PCR	PCR	
QM027-01	H12	PCR	PCR	PCR	