

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM029-01	A01	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM029-01	A02	MmiRQP0004	MIMAT0000522	mmu-let-7b-5p	
QM029-01	A03	MmiRQP0018	MIMAT0000655	mmu-miR-100-5p	
QM029-01	A04	MmiRQP0021	MIMAT0000133	mmu-miR-101a-3p	
QM029-01	A05	MmiRQP0032	MIMAT0000648	mmu-miR-10a-5p	
QM029-01	A06	MmiRQP0034	MIMAT0000208	mmu-miR-10b-5p	
QM029-01	A07	MmiRQP0056	MIMAT0000246	mmu-miR-122-5p	
QM029-01	A08	MmiRQP0074	MIMAT0000134	mmu-miR-124-3p	
QM029-01	A09	MmiRQP0096	MIMAT0000136	mmu-miR-125b-5p	
QM029-01	A10	MmiRQP0099	MIMAT0000138	mmu-miR-126-3p	
QM029-01	A11	MmiRQP0125	MIMAT0000140	mmu-miR-128-3p	
QM029-01	A12	MmiRQP0161	MIMAT0000144	mmu-miR-132-3p	
QM029-01	B01	MmiRQP0166	MIMAT0000145	mmu-miR-133a-3p	
QM029-01	B02	MmiRQP0167	MIMAT0000769	mmu-miR-133b-3p	
QM029-01	B03	MmiRQP0175	MIMAT0000149	mmu-miR-137-3p	
QM029-01	B04	MmiRQP0180	MIMAT0000656	mmu-miR-139-5p	
QM029-01	B05	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM029-01	B06	MmiRQP0184	MIMAT0000153	mmu-miR-141-3p	
QM029-01	B07	MmiRQP0186	MIMAT0000155	mmu-miR-142-3p	
QM029-01	B08	MmiRQP0188	MIMAT0000247	mmu-miR-143-3p	
QM029-01	B09	MmiRQP0192	MIMAT0000157	mmu-miR-145-5p	
QM029-01	B10	MmiRQP0196	MIMAT0000158	mmu-miR-146a-5p	
QM029-01	B11	MmiRQP0197	MIMAT0003475	mmu-miR-146b-5p	
QM029-01	B12	MmiRQP0204	MIMAT0000516	mmu-miR-148a-3p	
QM029-01	C01	MmiRQP0206	MIMAT0000580	mmu-miR-148b-3p	
QM029-01	C02	MmiRQP0212	MIMAT0004536	mmu-miR-151-5p	
QM029-01	C03	MmiRQP0223	MIMAT0000526	mmu-miR-15a-5p	
QM029-01	C04	MmiRQP0227	MIMAT0000527	mmu-miR-16-5p	
QM029-01	C05	MmiRQP0234	MIMAT0000673	mmu-miR-181b-5p	
QM029-01	C06	MmiRQP0245	MIMAT0000213	mmu-miR-184-3p	
QM029-01	C07	MmiRQP0248	MIMAT0000215	mmu-miR-186-5p	
QM029-01	C08	MmiRQP0255	MIMAT0000528	mmu-miR-18a-5p	
QM029-01	C09	MmiRQP0263	MIMAT0000221	mmu-miR-191-5p	
QM029-01	C10	MmiRQP0274	MIMAT0000517	mmu-miR-192-5p	
QM029-01	C11	MmiRQP0276	MIMAT0000223	mmu-miR-193-3p	
QM029-01	C12	MmiRQP0280	MIMAT0000224	mmu-miR-194-5p	
QM029-01	D01	MmiRQP0283	MIMAT0000225	mmu-miR-195-5p	
QM029-01	D02	MmiRQP0284	MIMAT0000518	mmu-miR-196a-5p	
QM029-01	D03	MmiRQP0290	MIMAT0000229	mmu-miR-199a-5p	
QM029-01	D04	MmiRQP0293	MIMAT0000651	mmu-miR-19a-3p	
QM029-01	D05	MmiRQP0044	MIMAT0000123	mmu-miR-1a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM029-01	D06	MmiRQP0298	MIMAT0000519	mmu-miR-200a-3p	
QM029-01	D07	MmiRQP0300	MIMAT0000233	mmu-miR-200b-3p	
QM029-01	D08	MmiRQP0305	MIMAT0000236	mmu-miR-203-3p	
QM029-01	D09	MmiRQP0306	MIMAT0000237	mmu-miR-204-5p	
QM029-01	D10	MmiRQP0307	MIMAT0000238	mmu-miR-205-5p	
QM029-01	D11	MmiRQP0317	MIMAT0000658	mmu-miR-210-3p	
QM029-01	D12	MmiRQP0320	MIMAT0000661	mmu-miR-214-3p	
QM029-01	E01	MmiRQP0324	MIMAT0003729	mmu-miR-216b-5p	
QM029-01	E02	MmiRQP0338	MIMAT0000669	mmu-miR-221-3p	
QM029-01	E03	MmiRQP0339	MIMAT0000670	mmu-miR-222-3p	
QM029-01	E04	MmiRQP0342	MIMAT0000665	mmu-miR-223-3p	
QM029-01	E05	MmiRQP0332	MIMAT0000531	mmu-miR-22-3p	
QM029-01	E06	MmiRQP0344	MIMAT0000532	mmu-miR-23a-3p	
QM029-01	E07	MmiRQP0346	MIMAT0000125	mmu-miR-23b-3p	
QM029-01	E08	MmiRQP0354	MIMAT0000533	mmu-miR-26a-5p	
QM029-01	E09	MmiRQP0357	MIMAT0000534	mmu-miR-26b-5p	
QM029-01	E10	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM029-01	E11	MmiRQP0361	MIMAT0000126	mmu-miR-27b-3p	
QM029-01	E12	MmiRQP0365	MIMAT0004576	mmu-miR-296-3p	
QM029-01	F01	MmiRQP0364	MIMAT0000374	mmu-miR-296-5p	
QM029-01	F02	MmiRQP0371	MIMAT0000535	mmu-miR-29a-3p	
QM029-01	F03	MmiRQP0373	MIMAT0000127	mmu-miR-29b-3p	
QM029-01	F04	MmiRQP0375	MIMAT0000536	mmu-miR-29c-3p	
QM029-01	F05	MmiRQP0381	MIMAT0000380	mmu-miR-302a-3p	
QM029-01	F06	MmiRQP0393	MIMAT0000130	mmu-miR-30b-5p	
QM029-01	F07	MmiRQP1224	MIMAT0000538	mmu-miR-31-5p	
QM029-01	F08	MmiRQP0418	MIMAT0004642	mmu-miR-330-5p	
QM029-01	F09	MmiRQP0420	MIMAT0000571	mmu-miR-331-3p	
QM029-01	F10	MmiRQP0419	MIMAT0004643	mmu-miR-331-5p	
QM029-01	F11	MmiRQP0426	MIMAT0000582	mmu-miR-338-3p	
QM029-01	F12	MmiRQP0434	MIMAT0004651	mmu-miR-340-5p	
QM029-01	G01	MmiRQP0436	MIMAT0000590	mmu-miR-342-3p	
QM029-01	G02	MmiRQP0440	MIMAT0000542	mmu-miR-34a-5p	
QM029-01	G03	MmiRQP0444	MIMAT0000381	mmu-miR-34c-5p	
QM029-01	G04	MmiRQP0456	MIMAT0001095	mmu-miR-370-3p	
QM029-01	G05	MmiRQP0464	MIMAT0003727	mmu-miR-374-5p	
QM029-01	G06	MmiRQP0466	MIMAT0000739	mmu-miR-375-3p	
QM029-01	G07	MmiRQP0473	MIMAT0003151	mmu-miR-378-3p	
QM029-01	G08	MmiRQP0536	MIMAT0003182	mmu-miR-494-3p	
QM029-01	G09	MmiRQP0786	MIMAT0000677	mmu-miR-7a-5p	
QM029-01	G10	MmiRQP0825	MIMAT0000142	mmu-miR-9-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM029-01	G11	MmiRQP0852	MIMAT0000541	mmu-miR-96-5p	
QM029-01	G12	MmiRQP0853	MIMAT0000545	mmu-miR-98-5p	
QM029-01	H01	NC	NC	NC	
QM029-01	H02	NC	NC	NC	
QM029-01	H03	MmiRQP9002	NR_003027	MK1	
QM029-01	H04	MmiRQP9012	NR_028128	MK2	
QM029-01	H05	MmiRQP9022	NR_028554	MK3	
QM029-01	H06	MmiRQP9042	NR_028550	MK4	
QM029-01	H07	MmiRQP9052	NR_028541	MK5	
QM029-01	H08	MmiRQP9062	NR_028091	MK6	
QM029-01	H09	RT	RT	RT	
QM029-01	H10	RT	RT	RT	
QM029-01	H11	PCR	PCR	PCR	
QM029-01	H12	PCR	PCR	PCR	