

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM020-01	A01	MmiRQP0096	MIMAT0000136	mmu-miR-125b-5p	
QM020-01	A02	MmiRQP0099	MIMAT0000138	mmu-miR-126-3p	
QM020-01	A03	MmiRQP0125	MIMAT0000140	mmu-miR-128-3p	
QM020-01	A04	MmiRQP0137	MIMAT0000209	mmu-miR-129-5p	
QM020-01	A05	MmiRQP0161	MIMAT0000144	mmu-miR-132-3p	
QM020-01	A06	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM020-01	A07	MmiRQP0192	MIMAT0000157	mmu-miR-145-5p	
QM020-01	A08	MmiRQP0208	MIMAT0000159	mmu-miR-149-5p	
QM020-01	A09	MmiRQP0223	MIMAT0000526	mmu-miR-15a-5p	
QM020-01	A10	MmiRQP0225	MIMAT0000124	mmu-miR-15b-5p	
QM020-01	A11	MmiRQP0227	MIMAT0000527	mmu-miR-16-5p	
QM020-01	A12	MmiRQP0248	MIMAT0000215	mmu-miR-186-5p	
QM020-01	B01	MmiRQP0283	MIMAT0000225	mmu-miR-195-5p	
QM020-01	B02	MmiRQP0290	MIMAT0000229	mmu-miR-199a-5p	
QM020-01	B03	MmiRQP0300	MIMAT0000233	mmu-miR-200b-3p	
QM020-01	B04	MmiRQP0302	MIMAT0000657	mmu-miR-200c-3p	
QM020-01	B05	MmiRQP0305	MIMAT0000236	mmu-miR-203-3p	
QM020-01	B06	MmiRQP0307	MIMAT0000238	mmu-miR-205-5p	
QM020-01	B07	MmiRQP0308	MIMAT0000239	mmu-miR-206-3p	
QM020-01	B08	MmiRQP0316	MIMAT0000530	mmu-miR-21-5p	
QM020-01	B09	MmiRQP0317	MIMAT0000658	mmu-miR-210-3p	
QM020-01	B10	MmiRQP0320	MIMAT0000661	mmu-miR-214-3p	
QM020-01	B11	MmiRQP0327	MIMAT0000663	mmu-miR-218-5p	
QM020-01	B12	MmiRQP0332	MIMAT0000531	mmu-miR-22-3p	
QM020-01	C01	MmiRQP0338	MIMAT0000669	mmu-miR-221-3p	
QM020-01	C02	MmiRQP0339	MIMAT0000670	mmu-miR-222-3p	
QM020-01	C03	MmiRQP0342	MIMAT0000665	mmu-miR-223-3p	
QM020-01	C04	MmiRQP0352	MIMAT0000652	mmu-miR-25-3p	
QM020-01	C05	MmiRQP0354	MIMAT0000533	mmu-miR-26a-5p	
QM020-01	C06	MmiRQP0357	MIMAT0000534	mmu-miR-26b-5p	
QM020-01	C07	MmiRQP0373	MIMAT0000127	mmu-miR-29b-3p	
QM020-01	C08	MmiRQP0415	MIMAT0000565	mmu-miR-328-3p	
QM020-01	C09	MmiRQP0434	MIMAT00004651	mmu-miR-340-5p	
QM020-01	C10	MmiRQP0520	MIMAT00003128	mmu-miR-485-5p	
QM020-01	C11	MmiRQP0537	MIMAT00003456	mmu-miR-495-3p	
QM020-01	C12	MmiRQP0837	MIMAT0000540	mmu-miR-93-5p	
QM020-01	D01	MmiRQP0852	MIMAT0000541	mmu-miR-96-5p	
QM020-01	D02	MmiRQP0890	MIMAT0000165	mmu-miR-155-5p	
QM020-01	D03	MmiRQP0895	MIMAT0000211	mmu-miR-182-5p	
QM020-01	D04	MmiRQP0918	MIMAT00004859	mmu-miR-193b-3p	
QM020-01	D05	MmiRQP0999	MIMAT00001537	mmu-miR-429-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM020-01	D06	MmiRQP1056	MIMAT0003112	mmu-miR-489-3p	
QM020-01	D07	MmiRQP1058	MIMAT0003453	mmu-miR-497-5p	
QM020-01	D08	MmiRQP1224	MIMAT0000538	mmu-miR-31-5p	
QM020-01	D09	MmiRQP0007	MIMAT0000383	mmu-let-7d-5p	
QM020-01	D10	MmiRQP0015	MIMAT0000121	mmu-let-7g-5p	
QM020-01	D11	MmiRQP0017	MIMAT0000122	mmu-let-7i-5p	
QM020-01	D12	MmiRQP0156	MIMAT0000141	mmu-miR-130a-3p	
QM020-01	E01	MmiRQP0159	MIMAT0000387	mmu-miR-130b-3p	
QM020-01	E02	MmiRQP0184	MIMAT0000153	mmu-miR-141-3p	
QM020-01	E03	MmiRQP0204	MIMAT0000516	mmu-miR-148a-3p	
QM020-01	E04	MmiRQP0255	MIMAT0000528	mmu-miR-18a-5p	
QM020-01	E05	MmiRQP0298	MIMAT0000519	mmu-miR-200a-3p	
QM020-01	E06	MmiRQP0312	MIMAT0000529	mmu-miR-20a-5p	
QM020-01	E07	MmiRQP0314	MIMAT0003187	mmu-miR-20b-5p	
QM020-01	E08	MmiRQP0853	MIMAT0000545	mmu-miR-98-5p	
QM020-01	E09	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM020-01	E10	MmiRQP0004	MIMAT0000522	mmu-let-7b-5p	
QM020-01	E11	MmiRQP0006	MIMAT0000523	mmu-let-7c-5p	
QM020-01	E12	MmiRQP0010	MIMAT0000524	mmu-let-7e-5p	
QM020-01	F01	MmiRQP0012	MIMAT0000525	mmu-let-7f-5p	
QM020-01	F02	MmiRQP0018	MIMAT0000655	mmu-miR-100-5p	
QM020-01	F03	MmiRQP0032	MIMAT0000648	mmu-miR-10a-5p	
QM020-01	F04	MmiRQP0034	MIMAT0000208	mmu-miR-10b-5p	
QM020-01	F05	MmiRQP0166	MIMAT0000145	mmu-miR-133a-3p	
QM020-01	F06	MmiRQP0230	MIMAT0000649	mmu-miR-17-5p	
QM020-01	F07	MmiRQP0232	MIMAT0000210	mmu-miR-181a-5p	
QM020-01	F08	MmiRQP0234	MIMAT0000673	mmu-miR-181b-5p	
QM020-01	F09	MmiRQP0235	MIMAT0000674	mmu-miR-181c-5p	
QM020-01	F10	MmiRQP0237	MIMAT0004324	mmu-miR-181d-5p	
QM020-01	F11	MmiRQP0293	MIMAT0000651	mmu-miR-19a-3p	
QM020-01	F12	MmiRQP0295	MIMAT0000513	mmu-miR-19b-3p	
QM020-01	G01	MmiRQP0306	MIMAT0000237	mmu-miR-204-5p	
QM020-01	G02	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM020-01	G03	MmiRQP0361	MIMAT0000126	mmu-miR-27b-3p	
QM020-01	G04	MmiRQP0371	MIMAT0000535	mmu-miR-29a-3p	
QM020-01	G05	MmiRQP0375	MIMAT0000536	mmu-miR-29c-3p	
QM020-01	G06	MmiRQP0786	MIMAT0000677	mmu-miR-7a-5p	
QM020-01	G07	MmiRQP0862	MIMAT0000385	mmu-miR-106a-5p	
QM020-01	G08	MmiRQP0289	MIMAT0004667	mmu-miR-199b-3p	mmu-miR-199a-3p
QM020-01	G09	MmiRQP0929	MIMAT0004628	mmu-miR-21-3p	
QM020-01	G10	MmiRQP0182	MIMAT0000152	mmu-miR-140-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM020-01	G11	MmiRQP0785	MIMAT0004670	mmu-miR-7a-1-3p	
QM020-01	G12	MmiRQP0921	MIMAT0004660	mmu-miR-19a-5p	
QM020-01	H01	NC	NC	NC	
QM020-01	H02	NC	NC	NC	
QM020-01	H03	MmiRQP9002	NR_003027	MK1	
QM020-01	H04	MmiRQP9012	NR_028128	MK2	
QM020-01	H05	MmiRQP9022	NR_028554	MK3	
QM020-01	H06	MmiRQP9042	NR_028550	MK4	
QM020-01	H07	MmiRQP9052	NR_028541	MK5	
QM020-01	H08	MmiRQP9062	NR_028091	MK6	
QM020-01	H09	RT	RT	RT	
QM020-01	H10	RT	RT	RT	
QM020-01	H11	PCR	PCR	PCR	
QM020-01	H12	PCR	PCR	PCR	