

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM021-01	A01	MmiRQP0002	MIMAT0000521	mmu-let-7a-5p	
QM021-01	A02	MmiRQP0004	MIMAT0000522	mmu-let-7b-5p	
QM021-01	A03	MmiRQP0006	MIMAT0000523	mmu-let-7c-5p	
QM021-01	A04	MmiRQP0029	MIMAT0000386	mmu-miR-106b-5p	
QM021-01	A05	MmiRQP0032	MIMAT0000648	mmu-miR-10a-5p	
QM021-01	A06	MmiRQP0074	MIMAT0000134	mmu-miR-124-3p	
QM021-01	A07	MmiRQP0093	MIMAT0004528	mmu-miR-125a-3p	
QM021-01	A08	MmiRQP0096	MIMAT0000136	mmu-miR-125b-5p	
QM021-01	A09	MmiRQP0099	MIMAT0000138	mmu-miR-126-3p	
QM021-01	A10	MmiRQP0111	MIMAT0000139	mmu-miR-127-3p	
QM021-01	A11	MmiRQP0125	MIMAT0000140	mmu-miR-128-3p	
QM021-01	A12	MmiRQP0139	MIMAT0000544	mmu-miR-129-2-3p	
QM021-01	B01	MmiRQP0137	MIMAT0000209	mmu-miR-129-5p	
QM021-01	B02	MmiRQP0159	MIMAT0000387	mmu-miR-130b-3p	
QM021-01	B03	MmiRQP0161	MIMAT0000144	mmu-miR-132-3p	
QM021-01	B04	MmiRQP0167	MIMAT0000769	mmu-miR-133b-3p	
QM021-01	B05	MmiRQP0172	MIMAT0000612	mmu-miR-135b-5p	
QM021-01	B06	MmiRQP0175	MIMAT0000149	mmu-miR-137-3p	
QM021-01	B07	MmiRQP0180	MIMAT0000656	mmu-miR-139-5p	
QM021-01	B08	MmiRQP0181	MIMAT0000151	mmu-miR-140-5p	
QM021-01	B09	MmiRQP0184	MIMAT0000153	mmu-miR-141-3p	
QM021-01	B10	MmiRQP0188	MIMAT0000247	mmu-miR-143-3p	
QM021-01	B11	MmiRQP0192	MIMAT0000157	mmu-miR-145-5p	
QM021-01	B12	MmiRQP0196	MIMAT0000158	mmu-miR-146a-5p	
QM021-01	C01	MmiRQP0210	MIMAT0000160	mmu-miR-150-5p	
QM021-01	C02	MmiRQP0225	MIMAT0000124	mmu-miR-15b-5p	
QM021-01	C03	MmiRQP0230	MIMAT0000649	mmu-miR-17-5p	
QM021-01	C04	MmiRQP0232	MIMAT0000210	mmu-miR-181a-5p	
QM021-01	C05	MmiRQP0244	MIMAT0000212	mmu-miR-183-5p	
QM021-01	C06	MmiRQP0255	MIMAT0000528	mmu-miR-18a-5p	
QM021-01	C07	MmiRQP0258	MIMAT0000220	mmu-miR-190-5p	
QM021-01	C08	MmiRQP0263	MIMAT0000221	mmu-miR-191-5p	
QM021-01	C09	MmiRQP0274	MIMAT0000517	mmu-miR-192-5p	
QM021-01	C10	MmiRQP0280	MIMAT0000224	mmu-miR-194-5p	
QM021-01	C11	MmiRQP0283	MIMAT0000225	mmu-miR-195-5p	
QM021-01	C12	MmiRQP0284	MIMAT0000518	mmu-miR-196a-5p	
QM021-01	D01	MmiRQP0290	MIMAT0000229	mmu-miR-199a-5p	
QM021-01	D02	MmiRQP0289	MIMAT0004667	mmu-miR-199b-3p	mmu-miR-199a-3p
QM021-01	D03	MmiRQP0293	MIMAT0000651	mmu-miR-19a-3p	
QM021-01	D04	MmiRQP0295	MIMAT0000513	mmu-miR-19b-3p	
QM021-01	D05	MmiRQP0044	MIMAT0000123	mmu-miR-1a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM021-01	D06	MmiRQP0298	MIMAT0000519	mmu-miR-200a-3p	
QM021-01	D07	MmiRQP0302	MIMAT0000657	mmu-miR-200c-3p	
QM021-01	D08	MmiRQP0305	MIMAT0000236	mmu-miR-203-3p	
QM021-01	D09	MmiRQP0306	MIMAT0000237	mmu-miR-204-5p	
QM021-01	D10	MmiRQP0312	MIMAT0000529	mmu-miR-20a-5p	
QM021-01	D11	MmiRQP0320	MIMAT0000661	mmu-miR-214-3p	
QM021-01	D12	MmiRQP0316	MIMAT0000530	mmu-miR-21-5p	
QM021-01	E01	MmiRQP0327	MIMAT0000663	mmu-miR-218-5p	
QM021-01	E02	MmiRQP0338	MIMAT0000669	mmu-miR-221-3p	
QM021-01	E03	MmiRQP0342	MIMAT0000665	mmu-miR-223-3p	
QM021-01	E04	MmiRQP0344	MIMAT0000532	mmu-miR-23a-3p	
QM021-01	E05	MmiRQP0346	MIMAT0000125	mmu-miR-23b-3p	
QM021-01	E06	MmiRQP0349	MIMAT0000219	mmu-miR-24-3p	
QM021-01	E07	MmiRQP0352	MIMAT0000652	mmu-miR-25-3p	
QM021-01	E08	MmiRQP0354	MIMAT0000533	mmu-miR-26a-5p	
QM021-01	E09	MmiRQP0359	MIMAT0000537	mmu-miR-27a-3p	
QM021-01	E10	MmiRQP0365	MIMAT0004576	mmu-miR-296-3p	
QM021-01	E11	MmiRQP0364	MIMAT0000374	mmu-miR-296-5p	
QM021-01	E12	MmiRQP0371	MIMAT0000535	mmu-miR-29a-3p	
QM021-01	F01	MmiRQP0373	MIMAT0000127	mmu-miR-29b-3p	
QM021-01	F02	MmiRQP0375	MIMAT0000536	mmu-miR-29c-3p	
QM021-01	F03	MmiRQP0391	MIMAT0000128	mmu-miR-30a-5p	
QM021-01	F04	MmiRQP0396	MIMAT0000514	mmu-miR-30c-5p	
QM021-01	F05	MmiRQP0400	MIMAT0000248	mmu-miR-30e-5p	
QM021-01	F06	MmiRQP1224	MIMAT0000538	mmu-miR-31-5p	
QM021-01	F07	MmiRQP0404	MIMAT0000654	mmu-miR-32-5p	
QM021-01	F08	MmiRQP0415	MIMAT0000565	mmu-miR-328-3p	
QM021-01	F09	MmiRQP0430	MIMAT0000667	mmu-miR-33-5p	
QM021-01	F10	MmiRQP0427	MIMAT0000584	mmu-miR-339-5p	
QM021-01	F11	MmiRQP0436	MIMAT0000590	mmu-miR-342-3p	
QM021-01	F12	MmiRQP0440	MIMAT0000542	mmu-miR-34a-5p	
QM021-01	G01	MmiRQP0444	MIMAT0000381	mmu-miR-34c-5p	
QM021-01	G02	MmiRQP0456	MIMAT0001095	mmu-miR-370-3p	
QM021-01	G03	MmiRQP0466	MIMAT0000739	mmu-miR-375-3p	
QM021-01	G04	MmiRQP0473	MIMAT0003151	mmu-miR-378-3p	
QM021-01	G05	MmiRQP0476	MIMAT0000743	mmu-miR-379-5p	
QM021-01	G06	MmiRQP0509	MIMAT0001632	mmu-miR-451	
QM021-01	G07	MmiRQP0786	MIMAT0000677	mmu-miR-7a-5p	
QM021-01	G08	MmiRQP0837	MIMAT0000540	mmu-miR-93-5p	
QM021-01	G09	MmiRQP0825	MIMAT0000142	mmu-miR-9-5p	
QM021-01	G10	MmiRQP0852	MIMAT0000541	mmu-miR-96-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM021-01	G11	MmiRQP0853	MIMAT0000545	mmu-miR-98-5p	
QM021-01	G12	MmiRQP0854	MIMAT0000131	mmu-miR-99a-5p	
QM021-01	H01	NC	NC	NC	
QM021-01	H02	NC	NC	NC	
QM021-01	H03	MmiRQP9002	NR_003027	MK1	
QM021-01	H04	MmiRQP9012	NR_028128	MK2	
QM021-01	H05	MmiRQP9022	NR_028554	MK3	
QM021-01	H06	MmiRQP9042	NR_028550	MK4	
QM021-01	H07	MmiRQP9052	NR_028541	MK5	
QM021-01	H08	MmiRQP9062	NR_028091	MK6	
QM021-01	H09	RT	RT	RT	
QM021-01	H10	RT	RT	RT	
QM021-01	H11	PCR	PCR	PCR	
QM021-01	H12	PCR	PCR	PCR	