

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM003-01	A01	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM003-01	A02	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM003-01	A03	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM003-01	A04	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM003-01	A05	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM003-01	A06	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM003-01	A07	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM003-01	A08	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM003-01	A09	HmiRQP0473	MIMAT0000732	hsa-miR-378a-3p	
QM003-01	A10	HmiRQP0600	MIMAT0002855	hsa-miR-520d-5p	
QM003-01	A11	HmiRQP0237	MIMAT0002821	hsa-miR-181d	
QM003-01	A12	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM003-01	B01	HmiRQP0139	MIMAT0004605	hsa-miR-129-2-3p	
QM003-01	B02	HmiRQP0109	MIMAT0005923	hsa-miR-1269a	
QM003-01	B03	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM003-01	B04	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM003-01	B05	HmiRQP0006	MIMAT0000064	hsa-let-7c	
QM003-01	B06	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM003-01	B07	HmiRQP0170	MIMAT0000428	hsa-miR-135a-5p	
QM003-01	B08	HmiRQP0578	MIMAT0002866	hsa-miR-517c-3p	
QM003-01	B09	HmiRQP0255	MIMAT0000072	hsa-miR-18a-5p	
QM003-01	B10	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM003-01	B11	HmiRQP0256	MIMAT0001412	hsa-miR-18b-5p	
QM003-01	B12	HmiRQP0598	MIMAT0002846	hsa-miR-520c-3p	
QM003-01	C01	HmiRQP0172	MIMAT0000758	hsa-miR-135b-5p	
QM003-01	C02	HmiRQP0597	MIMAT0002843	hsa-miR-520b	
QM003-01	C03	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM003-01	C04	HmiRQP0575	MIMAT0002852	hsa-miR-517a-3p	
QM003-01	C05	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM003-01	C06	HmiRQP0632	MIMAT0004812	hsa-miR-548d-5p	
QM003-01	C07	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM003-01	C08	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM003-01	C09	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM003-01	C10	HmiRQP0286	MIMAT0001080	hsa-miR-196b-5p	
QM003-01	C11	HmiRQP0601	MIMAT0002825	hsa-miR-520e	
QM003-01	C12	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM003-01	D01	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM003-01	D02	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM003-01	D03	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM003-01	D04	HmiRQP0638	MIMAT0005935	hsa-miR-548i	
QM003-01	D05	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM003-01	D06	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM003-01	D07	HmiRQP0630	MIMAT0004806	hsa-miR-548c-5p	
QM003-01	D08	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM003-01	D09	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM003-01	D10	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM003-01	D11	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM003-01	D12	HmiRQP0405	MIMAT0000510	hsa-miR-320a	
QM003-01	E01	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM003-01	E02	HmiRQP0166	MIMAT0000427	hsa-miR-133a	
QM003-01	E03	HmiRQP0591	MIMAT0005454	hsa-miR-519b-5p	hsa-miR-519c-5p
					hsa-miR-523-5p
					hsa-miR-518e-5p
					hsa-miR-522-5p
					hsa-miR-519a-5p
QM003-01	E04	HmiRQP0612	MIMAT0005455	hsa-miR-520c-5p	hsa-miR-526a
					hsa-miR-518d-5p
QM003-01	E05	HmiRQP0136	MIMAT0004548	hsa-miR-129-1-3p	
QM003-01	E06	HmiRQP0586	MIMAT0002841	hsa-miR-518f-5p	
QM003-01	E07	HmiRQP0390	MIMAT0000088	hsa-miR-30a-3p	
QM003-01	E08	HmiRQP2063	MIMAT0018954	hsa-miR-548ae	
QM003-01	E09	HmiRQP2098	MIMAT0018990	hsa-miR-548aj-3p	
QM003-01	E10	HmiRQP2166	MIMAT0019059	hsa-miR-1269b	
QM003-01	E11	HmiRQP2182	MIMAT0019074	hsa-miR-378i	
QM003-01	E12	HmiRQP2264	MIMAT0019732	hsa-miR-4662a-3p	
QM003-01	F01	HmiRQP2268	MIMAT0019736	hsa-miR-4662b	
QM003-01	F02				
QM003-01	F03				
QM003-01	F04				
QM003-01	F05				
QM003-01	F06				
QM003-01	F07				
QM003-01	F08				
QM003-01	F09				
QM003-01	F10				
QM003-01	F11				
QM003-01	F12				
QM003-01	G01				
QM003-01	G02				
QM003-01	G03				
QM003-01	G04				
QM003-01	G05				

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM003-01	G06				
QM003-01	G07				
QM003-01	G08				
QM003-01	G09				
QM003-01	G10				
QM003-01	G11				
QM003-01	G12				
QM003-01	H01	NC	NC	NC	
QM003-01	H02	NC	NC	NC	
QM003-01	H03	HmiRQP9001	NR_002752	HK1	
QM003-01	H04	HmiRQP9011	NR_002750	HK2	
QM003-01	H05	HmiRQP9021	NR_002745	HK3	
QM003-01	H06	HmiRQP9051	NR_002746	HK4	
QM003-01	H07	HmiRQP9061	NR_002744	HK5	
QM003-01	H08	HmiRQP9071	NR_002450	HK6	
QM003-01	H09	RT	RT	RT	
QM003-01	H10	RT	RT	RT	
QM003-01	H11	PCR	PCR	PCR	
QM003-01	H12	PCR	PCR	PCR	