

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM056_01	A01	HmiRQP0034	MIMAT0000254	hsa-mir-10b-5p	
QM056_01	A02	HmiRQP0225	MIMAT0000417	hsa-mir-15b-5p	
QM056_01	A03	HmiRQP0406	MIMAT0005792	hsa-mir-320b	
QM056_01	A04	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM056_01	A05	HmiRQP0172	MIMAT0000758	hsa-miR-135b-5p	
QM056_01	A06	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM056_01	A07	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM056_01	A08	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM056_01	A09	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM056_01	A10	HmiRQP0516	MIMAT0003150	hsa-miR-455-5p	
QM056_01	A11	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM056_01	A12	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM056_01	B01	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM056_01	B02	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM056_01	B03	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM056_01	B04	HmiRQP0404	MIMAT0000090	hsa-miR-32-5p	
QM056_01	B05	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM056_01	B06	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM056_01	B07	HmiRQP0213	MIMAT0000438	hsa-miR-152-3p	
QM056_01	B08	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	
QM056_01	B09	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM056_01	B10	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM056_01	B11	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM056_01	B12	HmiRQP0849	MIMAT0004987	hsa-miR-944	
QM056_01	C01	HmiRQP0037	MIMAT0005825	hsa-miR-1180-3p	
QM056_01	C02	HmiRQP0228	MIMAT0004518	hsa-miR-16-2-3p	
QM056_01	C03	HmiRQP0297	MIMAT0001620	hsa-miR-200a-5p	
QM056_01	C04	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM056_01	C05	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM056_01	C06	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM056_01	C07	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM056_01	C08	HmiRQP0078	MIMAT0005898	hsa-miR-1246	
QM056_01	C09	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM056_01	C10	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM056_01	C11	HmiRQP0317	MIMAT0000267	hsa-miR-210-3p	
QM056_01	C12	HmiRQP0402	MIMAT0004504	hsa-miR-31-3p	
QM056_01	D01	HmiRQP0510	MIMAT0001635	hsa-miR-452-5p	
QM056_01	E01	NC	NC	NC	
QM056_01	E02	NC	NC	NC	
QM056_01	E03	HKG1	HKG1	HKG1	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM056_01	E04	HKG2	HKG2	HKG2	
QM056_01	E05	HKG3	HKG3	HKG3	
QM056_01	E06	HKG4	HKG4	HKG4	
QM056_01	E07	HKG5	HKG5	HKG5	
QM056_01	E08	HKG6	HKG6	HKG6	
QM056_01	E09	RT	RT	RT	
QM056_01	E10	RT	RT	RT	
QM056_01	E11	PCR	PCR	PCR	
QM056_01	E12	PCR	PCR	PCR	