

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM064-1	A01	HmiRQP0033	MIMAT0004556	hsa-miR-10b-3p	
QM064-1	A02	HmiRQP0286	MIMAT0001080	hsa-miR-196b-5p	
QM064-1	A03	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM064-1	A04	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM064-1	A05	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM064-1	A06	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM064-1	A07	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM064-1	A08	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM064-1	A09	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM064-1	A10	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM064-1	A11	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM064-1	A12	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM064-1	B01	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM064-1	B02	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM064-1	B03	HmiRQP0537	MIMAT0002817	hsa-miR-495-3p	
QM064-1	C01	NC	NC	NC	
QM064-1	C02	NC	NC	NC	
QM064-1	C03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM064-1	C04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM064-1	C05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM064-1	C06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM064-1	C07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM064-1	C08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM064-1	C09	RT	RT	RT	
QM064-1	C10	RT	RT	RT	
QM064-1	C11	PCR	PCR	PCR	
QM064-1	C12	PCR	PCR	PCR	