

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM071-1	A01	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM071-1	A02	HmiRQP0785	MIMAT0004553	hsa-miR-7-1-3p	
QM071-1	A03	HmiRQP0537	MIMAT0002817	hsa-miR-495-3p	
QM071-1	A04	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM071-1	A05	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM071-1	A06	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM071-1	A07	HmiRQP0171	MIMAT0004698	hsa-miR-135b-3p	
QM071-1	A08	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM071-1	A09	HmiRQP0397	MIMAT0004551	hsa-miR-30d-3p	
QM071-1	A10	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM071-1	A11	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM071-1	A12	HmiRQP0394	MIMAT0004674	hsa-miR-30c-1-3p	
QM071-1	B01	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM071-1	B02	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM071-1	B03	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM071-1	B04	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM071-1	B05	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM071-1	B06	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM071-1	B07	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM071-1	B08	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM071-1	B09	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM071-1	B10	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM071-1	B11	HmiRQP0299	MIMAT0004571	hsa-miR-200b-5p	
QM071-1	B12	HmiRQP0539	MIMAT0004768	hsa-miR-497-3p	
QM071-1	C01	HmiRQP0656	MIMAT0004793	hsa-miR-556-3p	
QM071-1	C02	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM071-1	C03	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM071-1	D01	NC	NC	NC	
QM071-1	D02	NC	NC	NC	
QM071-1	D03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM071-1	D04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM071-1	D05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM071-1	D06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM071-1	D07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM071-1	D08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM071-1	D09	RT	RT	RT	
QM071-1	D10	RT	RT	RT	
QM071-1	D11	PCR	PCR	PCR	
QM071-1	D12	PCR	PCR	PCR	