

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM069-1	A01	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM069-1	A02	HmiRQP0617	MIMAT0003163	hsa-miR-539-5p	
QM069-1	A03	HmiRQP0785	MIMAT0004553	hsa-miR-7-1-3p	
QM069-1	A04	HmiRQP1321	MIMAT0010357	hsa-miR-670-5p	
QM069-1	A05	HmiRQP0380	MIMAT0000683	hsa-miR-302a-5p	
QM069-1	A06	HmiRQP0480	MIMAT0000737	hsa-miR-382-5p	
QM069-1	A07	HmiRQP0158	MIMAT0004680	hsa-miR-130b-5p	
QM069-1	A08	HmiRQP0258	MIMAT0000458	hsa-miR-190a-5p	
QM069-1	A09	HmiRQP0466	MIMAT0000728	hsa-miR-375-3p	
QM069-1	A10	HmiRQP2069	MIMAT0018961	hsa-miR-4443	
QM069-1	A11	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM069-1	A12	HmiRQP0411	MIMAT0000762	hsa-miR-324-3p	
QM069-1	B01	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM069-1	B02	HmiRQP0175	MIMAT0000429	hsa-miR-137-3p	
QM069-1	B03	HmiRQP0493	MIMAT0004749	hsa-miR-424-3p	
QM069-1	B04	HmiRQP0591	MIMAT0002831	hsa-miR-519c-5p	
QM069-1	B05	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM069-1	B06	HmiRQP0392	MIMAT0004589	hsa-miR-30b-3p	
QM069-1	B07	HmiRQP2354	MIMAT0019824	hsa-miR-4715-5p	
QM069-1	B08	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM069-1	B09	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM069-1	B10	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM069-1	B11	HmiRQP0254	MIMAT0002891	hsa-miR-18a-3p	
QM069-1	B12	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM069-1	C01	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM069-1	C02	HmiRQP0294	MIMAT0004491	hsa-miR-19b-1-5p	
QM069-1	C03	HmiRQP0826	MIMAT0004970	hsa-miR-920	
QM069-1	C04	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM069-1	C05	HmiRQP0358	MIMAT0004501	hsa-miR-27a-5p	
QM069-1	C06	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM069-1	C07	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM069-1	C08	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM069-1	C09	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM069-1	C10	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM069-1	D01	NC	NC	NC	
QM069-1	D02	NC	NC	NC	
QM069-1	D03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM069-1	D04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM069-1	D05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM069-1	D06	HsnoRNA U47	NR_002746	HK4(SNORD47)	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM069-1	D07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM069-1	D08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM069-1	D09	RT	RT	RT	
QM069-1	D10	RT	RT	RT	
QM069-1	D11	PCR	PCR	PCR	
QM069-1	D12	PCR	PCR	PCR	