

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM057_01	A01	HmiRQP0339	MIMAT0000279	hsa-mir-222-3p	
QM057_01	A02	HmiRQP0341	MIMAT0004570	hsa-mir-223	
QM057_01	A03	HmiRQP2779	MIMAT0024599	hsa-mir-6126	
QM057_01	A04	HmiRQP0315	MIMAT0004494	hsa-mir-21	
QM057_01	A05	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM057_01	A06	HmiRQP3416	MIMAT0022727	hsa-miR-1307-5p	
QM057_01	A07	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM057_01	A08	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM057_01	A09	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM057_01	A10	HmiRQP0154	MIMAT0005951	hsa-miR-1307-3p	
QM057_01	A11	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM057_01	A12	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM057_01	B01	HmiRQP4732	MIMAT0037313	hsa-miR-375-5p	
QM057_01	B02	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM057_01	B03	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM057_01	B04	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM057_01	B05	HmiRQP0316	MIMAT0000076	hsa-mir-21-5p	
QM057_01	B06	HmiRQP0371	MIMAT0000086	hsa-mir-29a-3p	
QM057_01	C01	NC	NC	NC	
QM057_01	C02	NC	NC	NC	
QM057_01	C03	HKG1	HKG1	HKG1	
QM057_01	C04	HKG2	HKG2	HKG2	
QM057_01	C05	HKG3	HKG3	HKG3	
QM057_01	C06	HKG4	HKG4	HKG4	
QM057_01	C07	HKG5	HKG5	HKG5	
QM057_01	C08	HKG6	HKG6	HKG6	
QM057_01	C09	RT	RT	RT	
QM057_01	C10	RT	RT	RT	
QM057_01	C11	PCR	PCR	PCR	
QM057_01	C12	PCR	PCR	PCR	