

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM061_01	A01	HmiRQP0295	MIMAT0000074	hsa-mir-19b-3p	
QM061_01	A02	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM061_01	A03	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM061_01	A04	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM061_01	A05	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM061_01	A06	HmiRQP1314	MIMAT0009197	hsa-miR-205-3p	
QM061_01	A07	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM061_01	A08	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM061_01	A09	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM061_01	A10	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM061_01	B01	NC	NC	NC	
QM061_01	B02	NC	NC	NC	
QM061_01	B03	HKG1	HKG1	HKG1	
QM061_01	B04	HKG2	HKG2	HKG2	
QM061_01	B05	HKG3	HKG3	HKG3	
QM061_01	B06	HKG4	HKG4	HKG4	
QM061_01	B07	HKG5	HKG5	HKG5	
QM061_01	B08	HKG6	HKG6	HKG6	
QM061_01	B09	RT	RT	RT	
QM061_01	B10	RT	RT	RT	
QM061_01	B11	PCR	PCR	PCR	
QM061_01	B12	PCR	PCR	PCR	