

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM062_01	A01	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM062_01	A02	HmiRQP0347	MIMAT0004587	hsa-miR-23b-5p	
QM062_01	A03	HmiRQP0325	MIMAT0000274	hsa-miR-217-5p	
QM062_01	A04	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM062_01	A05	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM062_01	A06	HmiRQP0034	MIMAT0000254	hsa-mir-10b-5p	
QM062_01	A07	HmiRQP0154	MIMAT0005951	hsa-mir-1307	
QM062_01	A08	HmiRQP0161	MIMAT0000426	hsa-mir-132-3p	
QM062_01	A09	HmiRQP0247	MIMAT0000455	hsa-mir-185-5p	
QM062_01	A10	HmiRQP0283	MIMAT0000461	hsa-mir-195-5p	
QM062_01	A11	HmiRQP0295	MIMAT0000074	hsa-mir-19b-3p	
QM062_01	A12	HmiRQP0311	MIMAT0004493	hsa-mir-20a-3p	
QM062_01	B01	HmiRQP0315	MIMAT0004494	hsa-mir-21	
QM062_01	B02	HmiRQP0364	MIMAT0000690	hsa-mir-296-5p	
QM062_01	B03	HmiRQP0156	MIMAT0000425	hsa-mir-130a	
QM062_01	B04	HmiRQP0546	MIMAT0002872	hsa-mir-501	
QM062_01	B05	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM062_01	B06	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM062_01	B07	HmiRQP0078	MIMAT0005898	hsa-miR-1246	
QM062_01	B08	HmiRQP2236	MIMAT0019704	hsa-miR-4644	
QM062_01	C01	NC	NC	NC	
QM062_01	C02	NC	NC	NC	
QM062_01	C03	HKG1	HKG1	HKG1	
QM062_01	C04	HKG2	HKG2	HKG2	
QM062_01	C05	HKG3	HKG3	HKG3	
QM062_01	C06	HKG4	HKG4	HKG4	
QM062_01	C07	HKG5	HKG5	HKG5	
QM062_01	C08	HKG6	HKG6	HKG6	
QM062_01	C09	RT	RT	RT	
QM062_01	C10	RT	RT	RT	
QM062_01	C11	PCR	PCR	PCR	
QM062_01	C12	PCR	PCR	PCR	