

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM058_01	A01	HmiRQP0099	MIMAT0000445	hsa-mir-126-3p	
QM058_01	A02	HmiRQP0210	MIMAT0000451	hsa-mir-150-5p	
QM058_01	A03	HmiRQP0225	MIMAT0000417	hsa-mir-15b-5p	
QM058_01	A04	HmiRQP0229	MIMAT0000071	hsa-mir-17-3p	
QM058_01	A05	HmiRQP0234	MIMAT0000257	hsa-mir-181b-5p	
QM058_01	A06	HmiRQP0263	MIMAT0000440	hsa-mir-191-5p	
QM058_01	A07	HmiRQP0300	MIMAT0000318	hsa-mir-200b-3p	
QM058_01	A08	HmiRQP0302	MIMAT0000617	hsa-mir-200c-3p	
QM058_01	A09	HmiRQP0307	MIMAT0000266	hsa-mir-205-5p	
QM058_01	A10	HmiRQP0315	MIMAT0004494	hsa-mir-21	
QM058_01	A11	HmiRQP0316	MIMAT0000076	hsa-mir-21-5p	
QM058_01	A12	HmiRQP0354	MIMAT0000082	hsa-mir-26a-5p	
QM058_01	B01	HmiRQP0373	MIMAT0000100	hsa-mir-29b-3p	
QM058_01	B02	HmiRQP2084	MIMAT0018976	hsa-mir-4454	
QM058_01	B03	HmiRQP0837	MIMAT0000093	hsa-mir-93-5p	
QM058_01	C01	NC	NC	NC	
QM058_01	C02	NC	NC	NC	
QM058_01	C03	HKG1	HKG1	HKG1	
QM058_01	C04	HKG2	HKG2	HKG2	
QM058_01	C05	HKG3	HKG3	HKG3	
QM058_01	C06	HKG4	HKG4	HKG4	
QM058_01	C07	HKG5	HKG5	HKG5	
QM058_01	C08	HKG6	HKG6	HKG6	
QM058_01	C09	RT	RT	RT	
QM058_01	C10	RT	RT	RT	
QM058_01	C11	PCR	PCR	PCR	
QM058_01	C12	PCR	PCR	PCR	