

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM060_01	A01	HmiRQP0183	MIMAT0004598	hsa-mir-141	
QM060_01	A02	HmiRQP0301	MIMAT0004657	hsa-mir-200c	
QM060_01	A03	HmiRQP0339	MIMAT0000279	hsa-mir-222	
QM060_01	A04	HmiRQP4680	MIMAT0037306	hsa-miR-103a-1-5p	
QM060_01	A05	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM060_01	A06	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM060_01	A07	HmiRQP0434	MIMAT0004692	hsa-miR-340-5p	
QM060_01	A08	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM060_01	A09	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM060_01	A10	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM060_01	A11	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM060_01	A12	HmiRQP4732	MIMAT0037313	hsa-miR-375-5p	
QM060_01	B01	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM060_01	B02	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM060_01	B03	HmiRQP0023	MIMAT0007402	hsa-miR-103b	
QM060_01	B04	HmiRQP0358	MIMAT0004501	hsa-miR-27a-5p	
QM060_01	B05	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM060_01	B06	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM060_01	B07	HmiRQP0264	MIMAT0001618	hsa-miR-191-3p	
QM060_01	B08	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM060_01	B09	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM060_01	B10	HmiRQP2192	MIMAT0019337	hsa-miR-3960	
QM060_01	B11	HmiRQP2128	MIMAT0019022	hsa-miR-4488	
QM060_01	C01	NC	NC	NC	
QM060_01	C02	NC	NC	NC	
QM060_01	C03	HKG1	HKG1	HKG1	
QM060_01	C04	HKG2	HKG2	HKG2	
QM060_01	C05	HKG3	HKG3	HKG3	
QM060_01	C06	HKG4	HKG4	HKG4	
QM060_01	C07	HKG5	HKG5	HKG5	
QM060_01	C08	HKG6	HKG6	HKG6	
QM060_01	C09	RT	RT	RT	
QM060_01	C10	RT	RT	RT	
QM060_01	C11	PCR	PCR	PCR	
QM060_01	C12	PCR	PCR	PCR	