

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM055_01	A01	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM055_01	A02	HmiRQP3416	MIMAT0022727	hsa-miR-1307-5p	
QM055_01	A03	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM055_01	A04	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM055_01	A05	HmiRQP0809	MIMAT0004949	hsa-miR-877-5p	
QM055_01	A06	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM055_01	A07	HmiRQP0159	MIMAT0000691	hsa-miR-130b-3p	
QM055_01	A08	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM055_01	A09	HmiRQP0491	MIMAT0001340	hsa-miR-423-3p	
QM055_01	A10	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM055_01	A11	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM055_01	A12	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM055_01	B01	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM055_01	B02	HmiRQP2450	MIMAT0019919	hsa-miR-4767	
QM055_01	B03	HmiRQP0835	MIMAT0003218	hsa-miR-92b-3p	
QM055_01	B04	HmiRQP1496	MIMAT0015041	hsa-miR-1260b	
QM055_01	B05	HmiRQP0193	MIMAT0006789	hsa-miR-1468-5p	
QM055_01	B06	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM055_01	B07	HmiRQP0678	MIMAT0003242	hsa-miR-577	
QM055_01	B08	HmiRQP0134	MIMAT0005880	hsa-miR-1290	
QM055_01	B09	HmiRQP0233	MIMAT0004558	hsa-miR-181a-2-3p	
QM055_01	B10	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM055_01	B11	HmiRQP4536	MIMAT0030019	hsa-miR-7704	
QM055_01	C01	NC	NC	NC	
QM055_01	C02	NC	NC	NC	
QM055_01	C03	HKG1	HKG1	HKG1	
QM055_01	C04	HKG2	HKG2	HKG2	
QM055_01	C05	HKG3	HKG3	HKG3	
QM055_01	C06	HKG4	HKG4	HKG4	
QM055_01	C07	HKG5	HKG5	HKG5	
QM055_01	C08	HKG6	HKG6	HKG6	
QM055_01	C09	RT	RT	RT	
QM055_01	C10	RT	RT	RT	
QM055_01	C11	PCR	PCR	PCR	
QM055_01	C12	PCR	PCR	PCR	