

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM049_01	A01	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM049_01	A02	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM049_01	A03	HmiRQP0355	MIMAT0004681	hsa-miR-26a-2-3p	
QM049_01	A04	HmiRQP0812	MIMAT0004947	hsa-miR-885-5p	
QM049_01	A05	HmiRQP0115	MIMAT0005926	hsa-miR-1273a	
QM049_01	A06	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	
QM049_01	A07	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM049_01	A08	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM049_01	A09	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM049_01	A10	HmiRQP0297	MIMAT0001620	hsa-miR-200a-5p	
QM049_01	A11	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM049_01	A12	HmiRQP2076	MIMAT0018967	hsa-miR-4448	
QM049_01	B01	HmiRQP0251	MIMAT0000262	hsa-miR-187-3p	
QM049_01	B02	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM049_01	B03	HmiRQP0404	MIMAT0000090	hsa-miR-32-5p	
QM049_01	B04	HmiRQP2159	MIMAT0019053	hsa-miR-4516	
QM049_01	B05	HmiRQP0276	MIMAT0000459	hsa-miR-193a-3p	
QM049_01	B06	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM049_01	B07	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM049_01	B08	HmiRQP0562	MIMAT0002808	hsa-miR-511-5p	
QM049_01	C01	NC	NC	NC	
QM049_01	C02	NC	NC	NC	
QM049_01	C03	HKG1	HKG1	HKG1	
QM049_01	C04	HKG2	HKG2	HKG2	
QM049_01	C05	HKG3	HKG3	HKG3	
QM049_01	C06	HKG4	HKG4	HKG4	
QM049_01	C07	HKG5	HKG5	HKG5	
QM049_01	C08	HKG6	HKG6	HKG6	
QM049_01	C09	RT	RT	RT	
QM049_01	C10	RT	RT	RT	
QM049_01	C11	PCR	PCR	PCR	
QM049_01	C12	PCR	PCR	PCR	