

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM052_01	A01	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM052_01	A02	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM052_01	A03	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM052_01	A04	HmiRQP0190	MIMAT0000436	hsa-miR-144-3p	
QM052_01	A05	HmiRQP0212	MIMAT0004697	hsa-miR-151a-5p	
QM052_01	A06	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM052_01	A07	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM052_01	A08	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM052_01	A09	HmiRQP0378	MIMAT0000688	hsa-miR-301a-3p	
QM052_01	A10	HmiRQP0193	MIMAT0006789	hsa-miR-1468-5p	
QM052_01	A11	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM052_01	A12	HmiRQP0540	MIMAT0002820	hsa-miR-497-5p	
QM052_01	B01	HmiRQP0198	MIMAT0004766	hsa-miR-146b-3p	
QM052_01	B02	HmiRQP0317	MIMAT0000267	hsa-miR-210-3p	
QM052_01	B03	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM052_01	B04	HmiRQP0134	MIMAT0005880	hsa-miR-1290	
QM052_01	B05	HmiRQP0233	MIMAT0004558	hsa-miR-181a-2-3p	
QM052_01	B06	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM052_01	B07	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM052_01	B08	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM052_01	B09	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM052_01	B10	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM052_01	B11	HmiRQP0399	MIMAT0000693	hsa-miR-30e-3p	
QM052_01	B12	HmiRQP0078	MIMAT0005898	hsa-miR-1246	
QM052_01	C01	HmiRQP0228	MIMAT0004518	hsa-miR-16-2-3p	
QM052_01	C02	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM052_01	C03	HmiRQP1496	MIMAT0015041	hsa-miR-1260b	
QM052_01	C04	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM052_01	C05	HmiRQP2128	MIMAT0019022	hsa-miR-4488	
QM052_01	D01	NC	NC	NC	
QM052_01	D02	NC	NC	NC	
QM052_01	D03	HKG1	HKG1	HKG1	
QM052_01	D04	HKG2	HKG2	HKG2	
QM052_01	D05	HKG3	HKG3	HKG3	
QM052_01	D06	HKG4	HKG4	HKG4	
QM052_01	D07	HKG5	HKG5	HKG5	
QM052_01	D08	HKG6	HKG6	HKG6	
QM052_01	D09	RT	RT	RT	
QM052_01	D10	RT	RT	RT	
QM052_01	D11	PCR	PCR	PCR	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM052_01	D12	PCR	PCR	PCR	