

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM053_01	A01	HmiRQP0276	MIMAT0000459	hsa-mir-193a-3p	
QM053_01	A02	HmiRQP0317	MIMAT0000267	hsa-mir-210-3p	
QM053_01	A03	HmiRQP0344	MIMAT0000078	hsa-mir-23a	
QM053_01	A04	HmiRQP0486	MIMAT0003329	hsa-mir-411	
QM053_01	A05	HmiRQP0486	MIMAT0003329	hsa-mir-411-5p	
QM053_01	A06	HmiRQP2626	MIMAT0022259	hsa-mir-5100	
QM053_01	A07	HmiRQP3407	MIMAT0022721	hsa-mir-1247-3p	
QM053_01	A08	HmiRQP0191	MIMAT0004601	hsa-mir-145	
QM053_01	A09	HmiRQP0315	MIMAT0004494	hsa-mir-21	
QM053_01	A10	HmiRQP0337	MIMAT0004568	hsa-mir-221	
QM053_01	A11	HmiRQP0339	MIMAT0000279	hsa-mir-222	
QM053_01	A12	HmiRQP0341	MIMAT0004570	hsa-mir-223	
QM053_01	B01	HmiRQP0344	MIMAT0000078	hsa-mir-23a-3p	
QM053_01	B02	HmiRQP0439	MIMAT0004557	hsa-mir-34a	
QM053_01	B03	HmiRQP0504	MIMAT0001541	hsa-mir-449a	
QM053_01	B04	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM053_01	B05	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM053_01	B06	HmiRQP0442	MIMAT0000685	hsa-miR-34b-5p	
QM053_01	B07	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM053_01	B08	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	
QM053_01	B09	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM053_01	B10	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM053_01	B11	HmiRQP0070	MIMAT0005591	hsa-miR-1236	
QM053_01	B12	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM053_01	C01	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM053_01	C02	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM053_01	C03	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM053_01	C04	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM053_01	C05	HmiRQP0097	MIMAT0004603	hsa-miR-125b-2-3p	
QM053_01	C06	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM053_01	C07	HmiRQP1295	MIMAT0009198	hsa-miR-224-3p	
QM053_01	D01	NC	NC	NC	
QM053_01	D02	NC	NC	NC	
QM053_01	D03	HKG1	HKG1	HKG1	
QM053_01	D04	HKG2	HKG2	HKG2	
QM053_01	D05	HKG3	HKG3	HKG3	
QM053_01	D06	HKG4	HKG4	HKG4	
QM053_01	D07	HKG5	HKG5	HKG5	
QM053_01	D08	HKG6	HKG6	HKG6	
QM053_01	D09	RT	RT	RT	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM053_01	D10	RT	RT	RT	
QM053_01	D11	PCR	PCR	PCR	
QM053_01	D12	PCR	PCR	PCR	