

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM059_01	A01	HmiRQP4732	MIMAT0037313	hsa-miR-375-5p	
QM059_01	A02	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM059_01	A03	HmiRQP0517	MIMAT0004761	hsa-miR-483-5p	
QM059_01	A04	HmiRQP0466	MIMAT0000728	hsa-miR-375	
QM059_01	A05	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM059_01	A06	HmiRQP0118	MIMAT0005929	hsa-miR-1275	
QM059_01	A07	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM059_01	A08	HmiRQP0134	MIMAT0005880	hsa-miR-1290	
QM059_01	A09	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM059_01	A10	HmiRQP0674	MIMAT0004795	hsa-miR-574-5p	
QM059_01	A11	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM059_01	A12	HmiRQP0028	MIMAT0004672	hsa-mir-106b	
QM059_01	B01	HmiRQP0229	MIMAT0000071	hsa-mir-17	
QM059_01	B02	HmiRQP0254	MIMAT0002891	hsa-mir-18A	
QM059_01	B03	HmiRQP0302	MIMAT0000617	hsa-mir-200c-3p	
QM059_01	B04	HmiRQP0311	MIMAT0004493	hsa-mir-20a	
QM059_01	B05	HmiRQP0836	MIMAT0004509	hsa-mir-93	
QM059_01	B06	HmiRQP0184	MIMAT0000432	hsa-mir-141-3p	
QM059_01	B07	HmiRQP0192	MIMAT0000437	hsa-mir-145-5p	
QM059_01	B08	HmiRQP0857	MIMAT0000689	hsa-mir-99b-5p	
QM059_01	B09	HmiRQP0058	MIMAT0005458	hsa-miR-1224-5p	
QM059_01	B10	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM059_01	C01	NC	NC	NC	
QM059_01	C02	NC	NC	NC	
QM059_01	C03	HKG1	HKG1	HKG1	
QM059_01	C04	HKG2	HKG2	HKG2	
QM059_01	C05	HKG3	HKG3	HKG3	
QM059_01	C06	HKG4	HKG4	HKG4	
QM059_01	C07	HKG5	HKG5	HKG5	
QM059_01	C08	HKG6	HKG6	HKG6	
QM059_01	C09	RT	RT	RT	
QM059_01	C10	RT	RT	RT	
QM059_01	C11	PCR	PCR	PCR	
QM059_01	C12	PCR	PCR	PCR	