

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM051_01	A01	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM051_01	A02	HmiRQP0486	MIMAT0003329	hsa-miR-411-5p	
QM051_01	A03	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM051_01	A04	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM051_01	A05	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM051_01	A06	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM051_01	A07	HmiRQP0392	MIMAT0004589	hsa-miR-30b-3p	
QM051_01	A08	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM051_01	A09	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM051_01	A10	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM051_01	A11	HmiRQP1574	MIMAT0016878	hsa-miR-4257	
QM051_01	A12	HmiRQP0394	MIMAT0004674	hsa-miR-30c-1-3p	
QM051_01	B01	HmiRQP0395	MIMAT0004550	hsa-miR-30c-2-3p	
QM051_01	B02	HmiRQP4680	MIMAT0037306	hsa-miR-103a-1-5p	
QM051_01	B03	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM051_01	B04	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM051_01	B05	HmiRQP0023	MIMAT0007402	hsa-miR-103b	
QM051_01	B06	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM051_01	B07	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM051_01	B08	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM051_01	B09	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM051_01	B10	HmiRQP0305	MIMAT0000264	hsa-miR-203a-3p	
QM051_01	B11	HmiRQP2343	MIMAT0019813	hsa-miR-203b-5p	
QM051_01	B12	HmiRQP2344	MIMAT0019814	hsa-miR-203b-3p	
QM051_01	C01	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM051_01	C02	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM051_01	C03	HmiRQP0340	MIMAT0004569	hsa-miR-222-5p	
QM051_01	C04	HmiRQP0313	MIMAT0004752	hsa-mir-20b	
QM051_01	C05	HmiRQP0399	MIMAT0000693	hsa-mir-30e-3p	
QM051_01	C06	HmiRQP0315	MIMAT0004494	hsa-mir-21	
QM051_01	C07	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM051_01	D01	NC	NC	NC	
QM051_01	D02	NC	NC	NC	
QM051_01	D03	HKG1	HKG1	HKG1	
QM051_01	D04	HKG2	HKG2	HKG2	
QM051_01	D05	HKG3	HKG3	HKG3	
QM051_01	D06	HKG4	HKG4	HKG4	
QM051_01	D07	HKG5	HKG5	HKG5	
QM051_01	D08	HKG6	HKG6	HKG6	
QM051_01	D09	RT	RT	RT	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM051_01	D10	RT	RT	RT	
QM051_01	D11	PCR	PCR	PCR	
QM051_01	D12	PCR	PCR	PCR	