

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM072-1	A01	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM072-1	A02	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM072-1	A03	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM072-1	A04	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM072-1	A05	HmiRQP0348	MIMAT0000079	hsa-miR-24-1-5p	
QM072-1	A06	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM072-1	A07	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM072-1	A08	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM072-1	A09	HmiRQP0166	MIMAT0000427	hsa-miR-133a-3p	
QM072-1	A10	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM072-1	A11	HmiRQP0403	MIMAT0004505	hsa-miR-32-3p	
QM072-1	A12	HmiRQP0831	MIMAT0004507	hsa-miR-92a-1-5p	
QM072-1	B01	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM072-1	B02	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM072-1	B03	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM072-1	B04	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM072-1	B05	HmiRQP0517	MIMAT0004761	hsa-miR-483-5p	
QM072-1	B06	HmiRQP0125	MIMAT0000424	hsa-miR-128-3p	
QM072-1	B07	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM072-1	B08	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM072-1	B09	HmiRQP0372	MIMAT0004514	hsa-miR-29b-1-5p	
QM072-1	B10	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM072-1	B11	HmiRQP0466	MIMAT0000728	hsa-miR-375-3p	
QM072-1	B12	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM072-1	C01	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM072-1	C02	HmiRQP0224	MIMAT0004586	hsa-miR-15b-3p	
QM072-1	C03	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM072-1	C04	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM072-1	C05	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM072-1	C06	HmiRQP0214	MIMAT0000439	hsa-miR-153-3p	
QM072-1	C07	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM072-1	C08	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM072-1	C09	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM072-1	C10	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM072-1	C11	HmiRQP0531	MIMAT0002807	hsa-miR-491-5p	
QM072-1	C12	HmiRQP0006	MIMAT0000064	hsa-let-7c-5p	
QM072-1	D01	HmiRQP0014	MIMAT0004584	hsa-let-7g-3p	
QM072-1	D02	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM072-1	D03	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM072-1	D04	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM072-1	D05	HmiRQP0563	MIMAT0002823	hsa-miR-512-3p	
QM072-1	D06	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM072-1	D07	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM072-1	E01	NC	NC	NC	
QM072-1	E02	NC	NC	NC	
QM072-1	E03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM072-1	E04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM072-1	E05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM072-1	E06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM072-1	E07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM072-1	E08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM072-1	E09	RT	RT	RT	
QM072-1	E10	RT	RT	RT	
QM072-1	E11	PCR	PCR	PCR	
QM072-1	E12	PCR	PCR	PCR	