

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM067-1	A01	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM067-1	A02	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM067-1	A03	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM067-1	A04	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM067-1	A05	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM067-1	A06	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM067-1	A07	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM067-1	A08	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM067-1	A09	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM067-1	A10	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM067-1	A11	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM067-1	A12	HmiRQP0224	MIMAT0004586	hsa-miR-15b-3p	
QM067-1	B01	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM067-1	B02	HmiRQP0169	MIMAT0004595	hsa-miR-135a-3p	
QM067-1	B03	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM067-1	B04	HmiRQP0360	MIMAT0004588	hsa-miR-27b-5p	
QM067-1	B05	HmiRQP3046	MIMAT0000250	hsa-miR-139-5p	
QM067-1	B06	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM067-1	B07	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM067-1	B08	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM067-1	B09	HmiRQP0009	MIMAT0004485	hsa-let-7e-3p	
QM067-1	B10	HmiRQP0044	MIMAT0000416	hsa-miR-1-3p	
QM067-1	B11	HmiRQP0125	MIMAT0000424	hsa-miR-128-3p	
QM067-1	B12	HmiRQP0449	MIMAT0000707	hsa-miR-363-3p	
QM067-1	C01	HmiRQP0539	MIMAT0004768	hsa-miR-497-3p	
QM067-1	C02	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM067-1	C03	HmiRQP0168	MIMAT0000447	hsa-miR-134-5p	
QM067-1	C04	HmiRQP0170	MIMAT0000428	hsa-miR-135a-5p	
QM067-1	C05	HmiRQP0480	MIMAT0000737	hsa-miR-382-5p	
QM067-1	C06	HmiRQP0318	MIMAT0000268	hsa-miR-211-5p	
QM067-1	C07	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM067-1	C08	HmiRQP0831	MIMAT0004507	hsa-miR-92a-1-5p	
QM067-1	C09	HmiRQP0098	MIMAT0000444	hsa-miR-126-5p	
QM067-1	C10	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM067-1	C11	HmiRQP0372	MIMAT0004514	hsa-miR-29b-1-5p	
QM067-1	C12	HmiRQP0254	MIMAT0002891	hsa-miR-18a-3p	
QM067-1	D01	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM067-1	D02	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM067-1	D03	HmiRQP0834	MIMAT0004792	hsa-miR-92b-5p	
QM067-1	D04	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM067-1	D05	HmiRQP0483	MIMAT0001638	hsa-miR-409-5p	
QM067-1	D06	HmiRQP0297	MIMAT0001620	hsa-miR-200a-5p	
QM067-1	D07	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM067-1	D08	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM067-1	D09	HmiRQP0154	MIMAT0005951	hsa-miR-1307-3p	
QM067-1	D10	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM067-1	D11	HmiRQP0486	MIMAT0003329	hsa-miR-411-5p	
QM067-1	D12	HmiRQP0262	MIMAT0004929	hsa-miR-190b-5p	
QM067-1	E01	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM067-1	E02	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM067-1	E03	HmiRQP0537	MIMAT0002817	hsa-miR-495-3p	
QM067-1	E04	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM067-1	E05	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM067-1	E06	HmiRQP0252	MIMAT0000457	hsa-miR-188-5p	
QM067-1	E07	HmiRQP0138	MIMAT0005943	hsa-miR-1292-5p	
QM067-1	E08	HmiRQP0562	MIMAT0002808	hsa-miR-511-5p	
QM067-1	E09	HmiRQP0425	MIMAT0004701	hsa-miR-338-5p	
QM067-1	E10	HmiRQP0481	MIMAT0000738	hsa-miR-383-5p	
QM067-1	E11	HmiRQP0437	MIMAT0000772	hsa-miR-345-5p	
QM067-1	E12	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM067-1	F01	HmiRQP0435	MIMAT0004694	hsa-miR-342-5p	
QM067-1	F02	HmiRQP0033	MIMAT0004556	hsa-miR-10b-3p	
QM067-1	F03	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM067-1	F04	HmiRQP0804	MIMAT0004911	hsa-miR-874-3p	
QM067-1	F05	HmiRQP0515	MIMAT0004784	hsa-miR-455-3p	
QM067-1	F06	HmiRQP1535	MIMAT0015085	hsa-miR-3200-3p	
QM067-1	F07	HmiRQP0213	MIMAT0000438	hsa-miR-152-3p	
QM067-1	F08	HmiRQP0519	MIMAT0002174	hsa-miR-484	
QM067-1	F09	HmiRQP0124	MIMAT0005939	hsa-miR-1281	
QM067-1	F10	HmiRQP0615	MIMAT0002888	hsa-miR-532-5p	
QM067-1	G01	NC	NC	NC	
QM067-1	G02	NC	NC	NC	
QM067-1	G03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM067-1	G04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM067-1	G05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM067-1	G06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM067-1	G07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM067-1	G08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM067-1	G09	RT	RT	RT	
QM067-1	G10	RT	RT	RT	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM067-1	G11	PCR	PCR	PCR	
QM067-1	G12	PCR	PCR	PCR	