

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM070-1	A01	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM070-1	A02	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM070-1	A03	HmiRQP0011	MIMAT0004486	hsa-let-7f-1-3p	
QM070-1	A04	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM070-1	A05	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM070-1	A06	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM070-1	A07	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM070-1	A08	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM070-1	A09	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM070-1	A10	HmiRQP0061	MIMAT0005576	hsa-miR-1226-5p	
QM070-1	A11	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM070-1	A12	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM070-1	B01	HmiRQP0158	MIMAT0004680	hsa-miR-130b-5p	
QM070-1	B02	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM070-1	B03	HmiRQP0175	MIMAT0000429	hsa-miR-137-3p	
QM070-1	B04	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM070-1	B05	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM070-1	B06	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM070-1	B07	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM070-1	B08	HmiRQP0213	MIMAT0000438	hsa-miR-152-3p	
QM070-1	B09	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM070-1	B10	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM070-1	B11	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM070-1	B12	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM070-1	C01	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM070-1	C02	HmiRQP0243	MIMAT0004560	hsa-miR-183-3p	
QM070-1	C03	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM070-1	C04	HmiRQP0252	MIMAT0000457	hsa-miR-188-5p	
QM070-1	C05	HmiRQP0254	MIMAT0002891	hsa-miR-18a-3p	
QM070-1	C06	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM070-1	C07	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM070-1	C08	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM070-1	C09	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM070-1	C10	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM070-1	C11	HmiRQP0294	MIMAT0004491	hsa-miR-19b-1-5p	
QM070-1	C12	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM070-1	D01	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM070-1	D02	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM070-1	D03	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM070-1	D04	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM070-1	D05	HmiRQP0317	MIMAT0000267	hsa-miR-210-3p	
QM070-1	D06	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM070-1	D07	HmiRQP0322	MIMAT0000272	hsa-miR-215-5p	
QM070-1	D08	HmiRQP0323	MIMAT0000273	hsa-miR-216a-5p	
QM070-1	D09	HmiRQP0324	MIMAT0004959	hsa-miR-216b-5p	
QM070-1	D10	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM070-1	D11	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM070-1	D12	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM070-1	E01	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM070-1	E02	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM070-1	E03	HmiRQP0348	MIMAT0000079	hsa-miR-24-1-5p	
QM070-1	E04	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM070-1	E05	HmiRQP0356	MIMAT0004500	hsa-miR-26b-3p	
QM070-1	E06	HmiRQP0372	MIMAT0004514	hsa-miR-29b-1-5p	
QM070-1	E07	HmiRQP0380	MIMAT0000683	hsa-miR-302a-5p	
QM070-1	E08	HmiRQP0390	MIMAT0000088	hsa-miR-30a-3p	
QM070-1	E09	HmiRQP0392	MIMAT0004589	hsa-miR-30b-3p	
QM070-1	E10	HmiRQP0394	MIMAT0004674	hsa-miR-30c-1-3p	
QM070-1	E11	HmiRQP0397	MIMAT0004551	hsa-miR-30d-3p	
QM070-1	E12	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM070-1	F01	HmiRQP0403	MIMAT0004505	hsa-miR-32-3p	
QM070-1	F02	HmiRQP0405	MIMAT0000510	hsa-miR-320a-3p	
QM070-1	F03	HmiRQP0417	MIMAT0000751	hsa-miR-330-3p	
QM070-1	F04	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM070-1	F05	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM070-1	F06	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM070-1	F07	HmiRQP0459	MIMAT0000724	hsa-miR-372-3p	
QM070-1	F08	HmiRQP0460	MIMAT0000725	hsa-miR-373-5p	
QM070-1	F09	HmiRQP0462	MIMAT0004688	hsa-miR-374a-3p	
QM070-1	F10	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM070-1	F11	HmiRQP0466	MIMAT0000728	hsa-miR-375-3p	
QM070-1	F12	HmiRQP0467	MIMAT0000729	hsa-miR-376a-3p	
QM070-1	G01	HmiRQP0469	MIMAT0002172	hsa-miR-376b-3p	
QM070-1	G02	HmiRQP0474	MIMAT0000731	hsa-miR-378a-5p	
QM070-1	G03	HmiRQP0489	MIMAT0003339	hsa-miR-421	
QM070-1	G04	HmiRQP0491	MIMAT0001340	hsa-miR-423-3p	
QM070-1	G05	HmiRQP0519	MIMAT0002174	hsa-miR-484	
QM070-1	G06	HmiRQP0548	MIMAT0004775	hsa-miR-502-3p	
QM070-1	G07	HmiRQP0550	MIMAT0002874	hsa-miR-503-5p	
QM070-1	G08	HmiRQP0554	MIMAT0002878	hsa-miR-506-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM070-1	G09	HmiRQP0563	MIMAT0002823	hsa-miR-512-3p	
QM070-1	G10	HmiRQP0570	MIMAT0002827	hsa-miR-515-3p	
QM070-1	G11	HmiRQP0573	MIMAT0002859	hsa-miR-516b-5p	
QM070-1	G12	HmiRQP0588	MIMAT0002869	hsa-miR-519a-3p	
QM070-1	H01	NC	NC	NC	
QM070-1	H02	NC	NC	NC	
QM070-1	H03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM070-1	H04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM070-1	H05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM070-1	H06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM070-1	H07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM070-1	H08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM070-1	H09	RT	RT	RT	
QM070-1	H10	RT	RT	RT	
QM070-1	H11	PCR	PCR	PCR	
QM070-1	H12	PCR	PCR	PCR	
QM070-2	A01	HmiRQP0719	MIMAT0004804	hsa-miR-615-5p	
QM070-2	A02	HmiRQP0740	MIMAT0003299	hsa-miR-630	
QM070-2	A03	HmiRQP0785	MIMAT0004553	hsa-miR-7-1-3p	
QM070-2	A04	HmiRQP0809	MIMAT0004949	hsa-miR-877-5p	
QM070-2	A05	HmiRQP0811	MIMAT0004948	hsa-miR-885-3p	
QM070-2	A06	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM070-2	A07	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM070-2	A08	HmiRQP0850	MIMAT0000094	hsa-miR-95-3p	
QM070-2	A09	HmiRQP0851	MIMAT0004510	hsa-miR-96-3p	
QM070-2	A10	HmiRQP0853	MIMAT0000096	hsa-miR-98-5p	
QM070-2	B01	NC	NC	NC	
QM070-2	B02	NC	NC	NC	
QM070-2	B03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM070-2	B04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM070-2	B05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM070-2	B06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM070-2	B07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM070-2	B08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM070-2	B09	RT	RT	RT	
QM070-2	B10	RT	RT	RT	
QM070-2	B11	PCR	PCR	PCR	
QM070-2	B12	PCR	PCR	PCR	