

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM074-1	A01	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM074-1	A02	HmiRQP0305	MIMAT0000264	hsa-miR-203a-3p	
QM074-1	A03	HmiRQP0425	MIMAT0004701	hsa-miR-338-5p	
QM074-1	A04	HmiRQP0024	MIMAT0004516	hsa-miR-105-3p	
QM074-1	A05	HmiRQP4584	MIMAT0030996	hsa-miR-8069	
QM074-1	A06	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM074-1	A07	HmiRQP0246	MIMAT0004611	hsa-miR-185-3p	
QM074-1	A08	HmiRQP2778	MIMAT0024598	hsa-miR-6125	
QM074-1	A09	HmiRQP2594	MIMAT0021043	hsa-miR-5010-5p	
QM074-1	A10	HmiRQP0098	MIMAT0000444	hsa-miR-126-5p	
QM074-1	A11	HmiRQP2322	MIMAT0019791	hsa-miR-4697-5p	
QM074-1	A12	HmiRQP2147	MIMAT0019041	hsa-miR-4505	
QM074-1	B01	HmiRQP2084	MIMAT0018976	hsa-miR-4454	
QM074-1	B02	HmiRQP0851	MIMAT0004510	hsa-miR-96-3p	
QM074-1	B03	HmiRQP0356	MIMAT0004500	hsa-miR-26b-3p	
QM074-1	B04	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM074-1	B05	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM074-1	B06	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM074-1	B07	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM074-1	B08	HmiRQP0224	MIMAT0004586	hsa-miR-15b-3p	
QM074-1	B09	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM074-1	B10	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM074-1	B11	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM074-1	B12	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM074-1	C01	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM074-1	C02	HmiRQP0539	MIMAT0004768	hsa-miR-497-3p	
QM074-1	C03	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM074-1	C04	HmiRQP0360	MIMAT0004588	hsa-miR-27b-5p	
QM074-1	C05	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM074-1	C06	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM074-1	C07	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM074-1	C08	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM074-1	C09	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM074-1	C10	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM074-1	C11	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM074-1	C12	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM074-1	D01	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM074-1	D02	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM074-1	D03	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM074-1	D04	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM074-1	D05	HmiRQP0171	MIMAT0004698	hsa-miR-135b-3p	
QM074-1	D06	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM074-1	D07	HmiRQP0500	MIMAT0002815	hsa-miR-432-3p	
QM074-1	D08	HmiRQP0475	MIMAT0004690	hsa-miR-379-3p	
QM074-1	D09	HmiRQP0218	MIMAT0000453	hsa-miR-154-3p	
QM074-1	D10	HmiRQP0462	MIMAT0004688	hsa-miR-374a-3p	
QM074-1	D11	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM074-1	D12	HmiRQP0166	MIMAT0000427	hsa-miR-133a-3p	
QM074-1	E01	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM074-1	E02	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM074-1	E03	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM074-1	E04	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM074-1	E05	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM074-1	E06	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM074-1	E07	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM074-1	E08	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM074-1	E09	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM074-1	E10	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM074-1	E11	HmiRQP0170	MIMAT0000428	hsa-miR-135a-5p	
QM074-1	E12	HmiRQP0318	MIMAT0000268	hsa-miR-211-5p	
QM074-1	F01	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM074-1	F02	HmiRQP0588	MIMAT0002869	hsa-miR-519a-3p	
QM074-1	F03	HmiRQP0467	MIMAT0000729	hsa-miR-376a-3p	
QM074-1	F04	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM074-1	F05	HmiRQP0719	MIMAT0004804	hsa-miR-615-5p	
QM074-1	F06	HmiRQP0780	MIMAT0003880	hsa-miR-671-5p	
QM074-1	F07	HmiRQP0466	MIMAT0000728	hsa-miR-375-3p	
QM074-1	F08	HmiRQP0658	MIMAT0003221	hsa-miR-557	
QM074-1	F09	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM074-1	F10	HmiRQP0414	MIMAT0000756	hsa-miR-326	
QM074-1	F11	HmiRQP0520	MIMAT0002175	hsa-miR-485-5p	
QM074-1	F12	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM074-1	G01	HmiRQP0489	MIMAT0003339	hsa-miR-421	
QM074-1	G02	HmiRQP0522	MIMAT0004762	hsa-miR-486-3p	
QM074-1	G03	HmiRQP0495	MIMAT0003393	hsa-miR-425-5p	
QM074-1	G04	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM074-1	G05	HmiRQP0445	MIMAT0004682	hsa-miR-361-3p	
QM074-1	G06	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM074-1	G07	HmiRQP0483	MIMAT0001638	hsa-miR-409-5p	
QM074-1	G08	HmiRQP0622	MIMAT0004954	hsa-miR-543	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM074-1	G09	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM074-1	G10	HmiRQP0536	MIMAT0002816	hsa-miR-494-3p	
QM074-1	G11	HmiRQP0803	MIMAT0004953	hsa-miR-873-5p	
QM074-1	G12	HmiRQP0542	MIMAT0004772	hsa-miR-499a-3p	
QM074-1	H01	NC	NC	NC	
QM074-1	H02	NC	NC	NC	
QM074-1	H03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM074-1	H04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM074-1	H05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM074-1	H06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM074-1	H07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM074-1	H08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM074-1	H09	RT	RT	RT	
QM074-1	H10	RT	RT	RT	
QM074-1	H11	PCR	PCR	PCR	
QM074-1	H12	PCR	PCR	PCR	
QM074-2	A01	HmiRQP0211	MIMAT0000757	hsa-miR-151a-3p	
QM074-2	A02	HmiRQP0683	MIMAT0004797	hsa-miR-582-3p	
QM074-2	A03	HmiRQP0125	MIMAT0000424	hsa-miR-128-3p	
QM074-2	A04	HmiRQP0617	MIMAT0003163	hsa-miR-539-5p	
QM074-2	A05	HmiRQP0362	MIMAT0000085	hsa-miR-28-5p	
QM074-2	A06	HmiRQP0703	MIMAT0003267	hsa-miR-599	
QM074-2	A07	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM074-2	A08	HmiRQP0734	MIMAT0003295	hsa-miR-626	
QM074-2	A09	HmiRQP0485	MIMAT0002171	hsa-miR-410-3p	
QM074-2	A10	HmiRQP0454	MIMAT0000721	hsa-miR-369-3p	
QM074-2	A11	HmiRQP0120	MIMAT0005933	hsa-miR-1277-3p	
QM074-2	A12	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM074-2	B01	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM074-2	B02	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM074-2	B03	HmiRQP0804	MIMAT0004911	hsa-miR-874-3p	
QM074-2	B04	HmiRQP0325	MIMAT0000274	hsa-miR-217-5p	
QM074-2	B05	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM074-2	B06	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM074-2	B07	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM074-2	B08	HmiRQP0323	MIMAT0000273	hsa-miR-216a-5p	
QM074-2	B09	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM074-2	B10	HmiRQP0276	MIMAT0000459	hsa-miR-193a-3p	
QM074-2	B11	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM074-2	B12	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM074-2	C01	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM074-2	C02	HmiRQP0177	MIMAT0000430	hsa-miR-138-5p	
QM074-2	C03	HmiRQP0173	MIMAT0000448	hsa-miR-136-5p	
QM074-2	C04	HmiRQP0168	MIMAT0000447	hsa-miR-134-5p	
QM074-2	C05	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM074-2	C06	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM074-2	C07	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM074-2	C08	HmiRQP0317	MIMAT0000267	hsa-miR-210-3p	
QM074-2	C09	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM074-2	C10	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM074-2	C11	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM074-2	C12	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM074-2	D01	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM074-2	D02	HmiRQP0330	MIMAT0000276	hsa-miR-219a-5p	
QM074-2	D03	HmiRQP0214	MIMAT0000439	hsa-miR-153-3p	
QM074-2	D04	HmiRQP0175	MIMAT0000429	hsa-miR-137-3p	
QM074-2	D05	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM074-2	D06	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM074-2	D07	HmiRQP0044	MIMAT0000416	hsa-miR-1-3p	
QM074-2	D08	HmiRQP0482	MIMAT0001075	hsa-miR-384	
QM074-2	D09	HmiRQP4576	MIMAT0030988	hsa-miR-8061	
QM074-2	E01	NC	NC	NC	
QM074-2	E02	NC	NC	NC	
QM074-2	E03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM074-2	E04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM074-2	E05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM074-2	E06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM074-2	E07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM074-2	E08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM074-2	E09	RT	RT	RT	
QM074-2	E10	RT	RT	RT	
QM074-2	E11	PCR	PCR	PCR	
QM074-2	E12	PCR	PCR	PCR	