

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM065-1	A01	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM065-1	A02	HmiRQP0305	MIMAT0000264	hsa-miR-203a-3p	
QM065-1	A03	HmiRQP3768	MIMAT0027572	hsa-miR-6780b-5p	
QM065-1	A04	HmiRQP0425	MIMAT0004701	hsa-miR-338-5p	
QM065-1	A05	HmiRQP0507	MIMAT0004910	hsa-miR-450b-3p	
QM065-1	A06	HmiRQP2777	MIMAT0024597	hsa-miR-6124	
QM065-1	A07	HmiRQP0711	MIMAT0003275	hsa-miR-607	
QM065-1	A08	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM065-1	A09	HmiRQP0014	MIMAT0004584	hsa-let-7g-3p	
QM065-1	A10	HmiRQP4541	MIMAT0030413	hsa-miR-4433b-5p	
QM065-1	A11	HmiRQP0246	MIMAT0004611	hsa-miR-185-3p	
QM065-1	A12	HmiRQP2775	MIMAT0023714	hsa-miR-6089	
QM065-1	B01	HmiRQP2776	MIMAT0023715	hsa-miR-6090	
QM065-1	B02	HmiRQP0059	MIMAT0005572	hsa-miR-1225-5p	
QM065-1	B03	HmiRQP0297	MIMAT0001620	hsa-miR-200a-5p	
QM065-1	B04	HmiRQP0254	MIMAT0002891	hsa-miR-18a-3p	
QM065-1	B05	HmiRQP1458	MIMAT0015002	hsa-miR-466	
QM065-1	B06	HmiRQP2630	MIMAT0022265	hsa-miR-548ar-5p	
QM065-1	B07	HmiRQP2618	MIMAT0021128	hsa-miR-5196-5p	
QM065-1	B08	HmiRQP0098	MIMAT0000444	hsa-miR-126-5p	
QM065-1	B09	HmiRQP2280	MIMAT0019749	hsa-miR-4669	
QM065-1	B10	HmiRQP2142	MIMAT0019036	hsa-miR-4500	
QM065-1	B11	HmiRQP2132	MIMAT0019026	hsa-miR-4491	
QM065-1	B12	HmiRQP2040	MIMAT0018932	hsa-miR-378f	
QM065-1	C01	HmiRQP2034	MIMAT0018925	hsa-miR-1268b	
QM065-1	C02	HmiRQP1560	MIMAT0016864	hsa-miR-4312	
QM065-1	C03	HmiRQP0526	MIMAT0002804	hsa-miR-488-5p	
QM065-1	C04	HmiRQP0358	MIMAT0004501	hsa-miR-27a-5p	
QM065-1	C05	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM065-1	C06	HmiRQP0313	MIMAT0004752	hsa-miR-20b-3p	
QM065-1	C07	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM065-1	C08	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM065-1	C09	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM065-1	C10	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM065-1	C11	HmiRQP0224	MIMAT0004586	hsa-miR-15b-3p	
QM065-1	C12	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM065-1	D01	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM065-1	D02	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM065-1	D03	HmiRQP0205	MIMAT0004699	hsa-miR-148b-5p	
QM065-1	D04	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM065-1	D05	HmiRQP0064	MIMAT0005582	hsa-miR-1228-5p	
QM065-1	D06	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM065-1	D07	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM065-1	D08	HmiRQP0785	MIMAT0004553	hsa-miR-7-1-3p	
QM065-1	D09	HmiRQP0544	MIMAT0002871	hsa-miR-500a-3p	
QM065-1	D10	HmiRQP0539	MIMAT0004768	hsa-miR-497-3p	
QM065-1	D11	HmiRQP0493	MIMAT0004749	hsa-miR-424-3p	
QM065-1	D12	HmiRQP0433	MIMAT0000750	hsa-miR-340-3p	
QM065-1	E01	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM065-1	E02	HmiRQP0360	MIMAT0004588	hsa-miR-27b-5p	
QM065-1	E03	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM065-1	E04	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM065-1	E05	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM065-1	E06	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM065-1	E07	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM065-1	E08	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM065-1	E09	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM065-1	E10	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM065-1	E11	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM065-1	E12	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM065-1	F01	HmiRQP0136	MIMAT0004548	hsa-miR-129-1-3p	
QM065-1	F02	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM065-1	F03	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM065-1	F04	HmiRQP0033	MIMAT0004556	hsa-miR-10b-3p	
QM065-1	F05	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM065-1	F06	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM065-1	F07	HmiRQP0299	MIMAT0004571	hsa-miR-200b-5p	
QM065-1	F08	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM065-1	F09	HmiRQP0171	MIMAT0004698	hsa-miR-135b-3p	
QM065-1	F10	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM065-1	F11	HmiRQP0471	MIMAT0004689	hsa-miR-377-5p	
QM065-1	F12	HmiRQP0738	MIMAT0003298	hsa-miR-629-3p	
QM065-1	G01	HmiRQP0462	MIMAT0004688	hsa-miR-374a-3p	
QM065-1	G02	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM065-1	G03	HmiRQP1323	MIMAT0010251	hsa-miR-449c-5p	
QM065-1	G04	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM065-1	G05	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM065-1	G06	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM065-1	G07	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM065-1	G08	HmiRQP0630	MIMAT0022740	hsa-miR-548am-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM065-1	G09	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM065-1	G10	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM065-1	G11	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM065-1	G12	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM065-1	C01	NC	NC	NC	
QM065-1	C02	NC	NC	NC	
QM065-1	C03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM065-1	C04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM065-1	C05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM065-1	C06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM065-1	C07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM065-1	C08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM065-1	C09	RT	RT	RT	
QM065-1	C10	RT	RT	RT	
QM065-1	C11	PCR	PCR	PCR	
QM065-1	C12	PCR	PCR	PCR	
QM065-2	A01	HmiRQP0640	MIMAT0005882	hsa-miR-548k	
QM065-2	A02	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM065-2	A03	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM065-2	A04	HmiRQP0170	MIMAT0000428	hsa-miR-135a-5p	
QM065-2	A05	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM065-2	A06	HmiRQP0467	MIMAT0000729	hsa-miR-376a-3p	
QM065-2	A07	HmiRQP0845	MIMAT0004983	hsa-miR-940	
QM065-2	A08	HmiRQP0780	MIMAT0003880	hsa-miR-671-5p	
QM065-2	A09	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM065-2	A10	HmiRQP0367	MIMAT0004901	hsa-miR-298	
QM065-2	A11	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM065-2	A12	HmiRQP0550	MIMAT0002874	hsa-miR-503-5p	
QM065-2	B01	HmiRQP0778	MIMAT0004952	hsa-miR-665	
QM065-2	B02	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM065-2	B03	HmiRQP0495	MIMAT0003393	hsa-miR-425-5p	
QM065-2	B04	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM065-2	B05	HmiRQP0131	MIMAT0005878	hsa-miR-1287-5p	
QM065-2	B06	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM065-2	B07	HmiRQP0483	MIMAT0001638	hsa-miR-409-5p	
QM065-2	B08	HmiRQP0751	MIMAT0003311	hsa-miR-641	
QM065-2	B09	HmiRQP0847	MIMAT0004985	hsa-miR-942-5p	
QM065-2	B10	HmiRQP0828	MIMAT0004972	hsa-miR-922	
QM065-2	B11	HmiRQP0728	MIMAT0003291	hsa-miR-622	
QM065-2	B12	HmiRQP0534	MIMAT0003161	hsa-miR-493-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM065-2	C01	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM065-2	C02	HmiRQP0536	MIMAT0002816	hsa-miR-494-3p	
QM065-2	C03	HmiRQP0615	MIMAT0002888	hsa-miR-532-5p	
QM065-2	C04	HmiRQP0303	MIMAT0002811	hsa-miR-202-3p	
QM065-2	C05	HmiRQP0452	MIMAT0000719	hsa-miR-367-3p	
QM065-2	C06	HmiRQP0435	MIMAT0004694	hsa-miR-342-5p	
QM065-2	C07	HmiRQP0676	MIMAT0004796	hsa-miR-576-3p	
QM065-2	C08	HmiRQP0118	MIMAT0005929	hsa-miR-1275	
QM065-2	C09	HmiRQP0542	MIMAT0004772	hsa-miR-499a-3p	
QM065-2	C10	HmiRQP0211	MIMAT0000757	hsa-miR-151a-3p	
QM065-2	C11	HmiRQP0538	MIMAT0002818	hsa-miR-496	
QM065-2	C12	HmiRQP0125	MIMAT0000424	hsa-miR-128-3p	
QM065-2	D01	HmiRQP0537	MIMAT0002817	hsa-miR-495-3p	
QM065-2	D02	HmiRQP0423	MIMAT0000754	hsa-miR-337-3p	
QM065-2	D03	HmiRQP0623	MIMAT0003164	hsa-miR-544a	
QM065-2	D04	HmiRQP0479	MIMAT0000736	hsa-miR-381-3p	
QM065-2	D05	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM065-2	D06	HmiRQP0761	MIMAT0003321	hsa-miR-651-5p	
QM065-2	D07	HmiRQP0652	MIMAT0003215	hsa-miR-552-3p	
QM065-2	D08	HmiRQP0362	MIMAT0000085	hsa-miR-28-5p	
QM065-2	D09	HmiRQP0703	MIMAT0003267	hsa-miR-599	
QM065-2	D10	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM065-2	D11	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM065-2	D12	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM065-2	E01	HmiRQP0804	MIMAT0004911	hsa-miR-874-3p	
QM065-2	E02	HmiRQP0531	MIMAT0002807	hsa-miR-491-5p	
QM065-2	E03	HmiRQP0706	MIMAT0003270	hsa-miR-602	
QM065-2	E04	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM065-2	E05	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM065-2	E06	HmiRQP0449	MIMAT0000707	hsa-miR-363-3p	
QM065-2	E07	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM065-2	E08	HmiRQP0325	MIMAT0000274	hsa-miR-217-5p	
QM065-2	E09	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM065-2	E10	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM065-2	E11	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM065-2	E12	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM065-2	F01	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM065-2	F02	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM065-2	F03	HmiRQP0252	MIMAT0000457	hsa-miR-188-5p	
QM065-2	F04	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM065-2	F05	HmiRQP0213	MIMAT0000438	hsa-miR-152-3p	
QM065-2	F06	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM065-2	F07	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM065-2	F08	HmiRQP0177	MIMAT0000430	hsa-miR-138-5p	
QM065-2	F09	HmiRQP0168	MIMAT0000447	hsa-miR-134-5p	
QM065-2	F10	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM065-2	F11	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM065-2	F12	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM065-2	G01	HmiRQP0317	MIMAT0000267	hsa-miR-210-3p	
QM065-2	G02	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM065-2	G03	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM065-2	G04	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM065-2	G05	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM065-2	G06	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM065-2	G07	HmiRQP0214	MIMAT0000439	hsa-miR-153-3p	
QM065-2	G08	HmiRQP3046	MIMAT0000250	hsa-miR-139-5p	
QM065-2	G09	HmiRQP0175	MIMAT0000429	hsa-miR-137-3p	
QM065-2	G10	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM065-2	G11	HmiRQP0057	MIMAT0005459	hsa-miR-1224-3p	
QM065-2	G12	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM065-2	C01	NC	NC	NC	
QM065-2	C02	NC	NC	NC	
QM065-2	C03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM065-2	C04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM065-2	C05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM065-2	C06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM065-2	C07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM065-2	C08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM065-2	C09	RT	RT	RT	
QM065-2	C10	RT	RT	RT	
QM065-2	C11	PCR	PCR	PCR	
QM065-2	C12	PCR	PCR	PCR	
QM065-3	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1-3p	
QM065-3	A02	HmiRQP1443	MIMAT0014985	hsa-miR-3123	
QM065-3	A03	HmiRQP0853	MIMAT0000096	hsa-miR-98-5p	
QM065-3	A04	HmiRQP3702	MIMAT0027506	hsa-miR-6803-5p	
QM065-3	A05	HmiRQP1615	MIMAT0016919	hsa-miR-4292	
QM065-3	A06	HmiRQP3830	MIMAT0027634	hsa-miR-6867-5p	
QM065-3	A07	HmiRQP3688	MIMAT0027492	hsa-miR-6796-5p	
QM065-3	B01	NC	NC	NC	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM065-3	B02	NC	NC	NC	
QM065-3	B03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM065-3	B04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM065-3	B05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM065-3	B06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM065-3	B07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM065-3	B08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM065-3	B09	RT	RT	RT	
QM065-3	B10	RT	RT	RT	
QM065-3	B11	PCR	PCR	PCR	
QM065-3	B12	PCR	PCR	PCR	