

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM073-1	A01	HmiRQP0195	MIMAT0004608	hsa-mir-146a	
QM073-1	A02	HmiRQP3577	MIMAT0027381	hsa-mir-6740	
QM073-1	A03	HmiRQP0425	MIMAT0004701	hsa-mir-338	
QM073-1	A04	HmiRQP3569	MIMAT0027373	hsa-mir-6736	
QM073-1	A05	HmiRQP3565	MIMAT0027369	hsa-mir-6734	
QM073-1	A06	HmiRQP3652	MIMAT0027456	hsa-mir-6778	
QM073-1	A07	HmiRQP3846	MIMAT0027650	hsa-mir-6875	
QM073-1	A08	HmiRQP2785	MIMAT0024615	hsa-mir-6131	
QM073-1	A09	HmiRQP2989	MIMAT0025844	hsa-mir-6716	
QM073-1	A10	HmiRQP1621	MIMAT0016925	hsa-mir-500b	
QM073-1	A11	HmiRQP0282	MIMAT0004615	hsa-mir-195	
QM073-1	A12	HmiRQP0014	MIMAT0004584	hsa-let-7g	
QM073-1	B01	HmiRQP3786	MIMAT0027590	hsa-mir-6845	
QM073-1	B02	HmiRQP0246	MIMAT0004611	hsa-mir-185	
QM073-1	B03	HmiRQP0297	MIMAT0001620	hsa-mir-200a	
QM073-1	B04	HmiRQP2636	MIMAT0022271	hsa-mir-664b	
QM073-1	B05	HmiRQP0460	MIMAT0000725	hsa-mir-373	
QM073-1	B06	HmiRQP0098	MIMAT0000444	hsa-mir-126	
QM073-1	B07	HmiRQP2487	MIMAT0019956	hsa-mir-4787	
QM073-1	B08	HmiRQP2451	MIMAT0019920	hsa-mir-4768	
QM073-1	B09	HmiRQP2413	MIMAT0019882	hsa-mir-4747	
QM073-1	B10	HmiRQP2369	MIMAT0019838	hsa-mir-4723	
QM073-1	B11	HmiRQP2366	MIMAT0019836	hsa-mir-4722	
QM073-1	B12	HmiRQP2250	MIMAT0019718	hsa-mir-4653	
QM073-1	C01	HmiRQP2138	MIMAT0019032	hsa-mir-4497	
QM073-1	C02	HmiRQP2133	MIMAT0019027	hsa-mir-4492	
QM073-1	C03	HmiRQP2127	MIMAT0019021	hsa-mir-4487	
QM073-1	C04	HmiRQP2077	MIMAT0018968	hsa-mir-4449	
QM073-1	C05	HmiRQP2060	MIMAT0018951	hsa-mir-4435-1	
QM073-1	C06	HmiRQP2043	MIMAT0018935	hsa-mir-4422	
QM073-1	C07	HmiRQP0851	MIMAT0004510	hsa-mir-96	
QM073-1	C08	HmiRQP0358	MIMAT0004501	hsa-mir-27a	
QM073-1	C09	HmiRQP0356	MIMAT0004500	hsa-mir-26b	
QM073-1	C10	HmiRQP0315	MIMAT0004494	hsa-mir-21	
QM073-1	C11	HmiRQP0313	MIMAT0004752	hsa-mir-20b	
QM073-1	C12	HmiRQP0238	MIMAT0000260	hsa-mir-182	
QM073-1	D01	HmiRQP0229	MIMAT0000071	hsa-mir-17	
QM073-1	D02	HmiRQP0224	MIMAT0004586	hsa-mir-15b	
QM073-1	D03	HmiRQP0222	MIMAT0004488	hsa-mir-15a	
QM073-1	D04	HmiRQP0209	MIMAT0004610	hsa-mir-150	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM073-1	D05	HmiRQP0207	MIMAT0004609	hsa-mir-149	
QM073-1	D06	HmiRQP0095	MIMAT0004592	hsa-mir-125b-1	
QM073-1	D07	HmiRQP1612	MIMAT0016916	hsa-mir-4286	
QM073-1	D08	HmiRQP1518	MIMAT0015066	hsa-mir-3065	
QM073-1	D09	HmiRQP1460	MIMAT0015004	hsa-mir-544b	
QM073-1	D10	HmiRQP0836	MIMAT0004509	hsa-mir-93	
QM073-1	D11	HmiRQP0824	MIMAT0000442	hsa-mir-9-1	
QM073-1	D12	HmiRQP0539	MIMAT0004768	hsa-mir-497	
QM073-1	E01	HmiRQP0513	MIMAT0003884	hsa-mir-454	
QM073-1	E02	HmiRQP0493	MIMAT0004749	hsa-mir-424	
QM073-1	E03	HmiRQP0474	MIMAT0000731	hsa-mir-378a	
QM073-1	E04	HmiRQP0433	MIMAT0000750	hsa-mir-340	
QM073-1	E05	HmiRQP0372	MIMAT0004514	hsa-mir-29b-1	
QM073-1	E06	HmiRQP0370	MIMAT0004503	hsa-mir-29a	
QM073-1	E07	HmiRQP0360	MIMAT0004588	hsa-mir-27b	
QM073-1	E08	HmiRQP0351	MIMAT0004498	hsa-mir-25	
QM073-1	E09	HmiRQP0341	MIMAT0004570	hsa-mir-223	
QM073-1	E10	HmiRQP0311	MIMAT0004493	hsa-mir-20a	
QM073-1	E11	HmiRQP0301	MIMAT0004657	hsa-mir-200c	
QM073-1	E12	HmiRQP0220	MIMAT0004658	hsa-mir-155	
QM073-1	F01	HmiRQP0203	MIMAT0004549	hsa-mir-148a	
QM073-1	F02	HmiRQP0189	MIMAT0004600	hsa-mir-144	
QM073-1	F03	HmiRQP0187	MIMAT0004599	hsa-mir-143	
QM073-1	F04	HmiRQP0160	MIMAT0004594	hsa-mir-132	
QM073-1	F05	HmiRQP0073	MIMAT0004591	hsa-mir-124-1	
QM073-1	F06	HmiRQP0055	MIMAT0004590	hsa-mir-122	
QM073-1	F07	HmiRQP0033	MIMAT0004556	hsa-mir-10b	
QM073-1	F08	HmiRQP0028	MIMAT0004672	hsa-mir-106b	
QM073-1	F09	HmiRQP0016	MIMAT0004585	hsa-let-7i	
QM073-1	F10	HmiRQP0009	MIMAT0004485	hsa-let-7e	
QM073-1	F11	HmiRQP0001	MIMAT0004481	hsa-let-7a-1	
QM073-1	F12	HmiRQP0299	MIMAT0004571	hsa-mir-200b	
QM073-1	G01	HmiRQP0439	MIMAT0004557	hsa-mir-34a	
QM073-1	G02	HmiRQP0498	MIMAT0004757	hsa-mir-431	
QM073-1	G03	HmiRQP0462	MIMAT0004688	hsa-mir-374a	
QM073-1	G04	HmiRQP0289	MIMAT0000232	hsa-mir-199a-1	
QM073-1	G05	HmiRQP1297	MIMAT0011158	hsa-mir-2115	
QM073-1	G06	HmiRQP0260	MIMAT0007883	hsa-mir-1909	
QM073-1	G07	HmiRQP0344	MIMAT0000078	hsa-mir-23a	
QM073-1	G08	HmiRQP0234	MIMAT0000257	hsa-mir-181b-1	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM073-1	G09	HmiRQP0295	MIMAT0000074	hsa-mir-19b-1	
QM073-1	G10	HmiRQP0167	MIMAT0000770	hsa-mir-133b	
QM073-1	G11	HmiRQP0630	MIMAT0022740	hsa-mir-548am	
QM073-1	G12	HmiRQP0030	MIMAT0000104	hsa-mir-107	
QM073-1	H01	NC	NC	NC	
QM073-1	H02	NC	NC	NC	
QM073-1	H03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM073-1	H04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM073-1	H05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM073-1	H06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM073-1	H07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM073-1	H08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM073-1	H09	RT	RT	RT	
QM073-1	H10	RT	RT	RT	
QM073-1	H11	PCR	PCR	PCR	
QM073-1	H12	PCR	PCR	PCR	
QM073-2	A01	HmiRQP0022	MIMAT0000101	hsa-mir-103a-1	
QM073-2	A02	HmiRQP0232	MIMAT0000256	hsa-mir-181a-1	
QM073-2	A03	HmiRQP0286	MIMAT0001080	hsa-mir-196b	
QM073-2	A04	HmiRQP0375	MIMAT0000681	hsa-mir-29c	
QM073-2	A05	HmiRQP0235	MIMAT0000258	hsa-mir-181c	
QM073-2	A06	HmiRQP0026	MIMAT0000103	hsa-mir-106a	
QM073-2	A07	HmiRQP0256	MIMAT0001412	hsa-mir-18b	
QM073-2	A08	HmiRQP0284	MIMAT0000226	hsa-mir-196a-1	
QM073-2	A09	HmiRQP0170	MIMAT0000428	hsa-mir-135a-1	
QM073-2	A10	HmiRQP0548	MIMAT0004775	hsa-mir-502	
QM073-2	A11	HmiRQP0318	MIMAT0000268	hsa-mir-211	
QM073-2	A12	HmiRQP0306	MIMAT0000265	hsa-mir-204	
QM073-2	B01	HmiRQP0592	MIMAT0002853	hsa-mir-519d	
QM073-2	B02	HmiRQP0274	MIMAT0000222	hsa-mir-192	
QM073-2	B03	HmiRQP0469	MIMAT0002172	hsa-mir-376b	
QM073-2	B04	HmiRQP0007	MIMAT0000065	hsa-let-7d	
QM073-2	B05	HmiRQP0464	MIMAT0004955	hsa-mir-374b	
QM073-2	B06	HmiRQP0772	MIMAT0003324	hsa-mir-661	
QM073-2	B07	HmiRQP0719	MIMAT0004804	hsa-mir-615	
QM073-2	B08	HmiRQP0438	MIMAT0000773	hsa-mir-346	
QM073-2	B09	HmiRQP0780	MIMAT0003880	hsa-mir-671	
QM073-2	B10	HmiRQP0466	MIMAT0000728	hsa-mir-375	
QM073-2	B11	HmiRQP0718	MIMAT0003282	hsa-mir-614	
QM073-2	B12	HmiRQP0367	MIMAT0004901	hsa-mir-298	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM073-2	C01	HmiRQP0287	MIMAT0000227	hsa-mir-197	
QM073-2	C02	HmiRQP0673	MIMAT0003239	hsa-mir-574	
QM073-2	C03	HmiRQP0584	MIMAT0002864	hsa-mir-518d	
QM073-2	C04	HmiRQP0760	MIMAT0003320	hsa-mir-650	
QM073-2	C05	HmiRQP0550	MIMAT0002874	hsa-mir-503	
QM073-2	C06	HmiRQP0068	MIMAT0005588	hsa-mir-1233-1	
QM073-2	C07	HmiRQP0066	MIMAT0005584	hsa-mir-1229	
QM073-2	C08	HmiRQP0414	MIMAT0000756	hsa-mir-326	
QM073-2	C09	HmiRQP0520	MIMAT0002175	hsa-mir-485	
QM073-2	C10	HmiRQP0811	MIMAT0004948	hsa-mir-885	
QM073-2	C11	HmiRQP0093	MIMAT0004602	hsa-mir-125a	
QM073-2	C12	HmiRQP0649	MIMAT0003214	hsa-mir-551a	
QM073-2	D01	HmiRQP0495	MIMAT0003393	hsa-mir-425	
QM073-2	D02	HmiRQP0443	MIMAT0004677	hsa-mir-34c	
QM073-2	D03	HmiRQP0502	MIMAT0001627	hsa-mir-433	
QM073-2	D04	HmiRQP0517	MIMAT0004761	hsa-mir-483	
QM073-2	D05	HmiRQP0445	MIMAT0004682	hsa-mir-361	
QM073-2	D06	HmiRQP0320	MIMAT0000271	hsa-mir-214	
QM073-2	D07	HmiRQP0113	MIMAT0005796	hsa-mir-1271	
QM073-2	D08	HmiRQP0809	MIMAT0004949	hsa-mir-877	
QM073-2	D09	HmiRQP0546	MIMAT0002872	hsa-mir-501	
QM073-2	D10	HmiRQP0756	MIMAT0003316	hsa-mir-646	
QM073-2	D11	HmiRQP0346	MIMAT0000418	hsa-mir-23b	
QM073-2	D12	HmiRQP0615	MIMAT0002888	hsa-mir-532	
QM073-2	E01	HmiRQP0303	MIMAT0002811	hsa-mir-202	
QM073-2	E02	HmiRQP0153	MIMAT0005950	hsa-mir-1306	
QM073-2	E03	HmiRQP0699	MIMAT0003263	hsa-mir-595	
QM073-2	E04	HmiRQP0717	MIMAT0003281	hsa-mir-613	
QM073-2	E05	HmiRQP0515	MIMAT0004784	hsa-mir-455	
QM073-2	E06	HmiRQP0551	MIMAT0002875	hsa-mir-504	
QM073-2	E07	HmiRQP0666	MIMAT0003231	hsa-mir-567	
QM073-2	E08	HmiRQP0078	MIMAT0005898	hsa-mir-1246	
QM073-2	E09	HmiRQP0435	MIMAT0004694	hsa-mir-342	
QM073-2	E10	HmiRQP0676	MIMAT0004796	hsa-mir-576	
QM073-2	E11	HmiRQP0481	MIMAT0000738	hsa-mir-383	
QM073-2	E12	HmiRQP0562	MIMAT0002808	hsa-mir-511	
QM073-2	F01	HmiRQP0211	MIMAT0000757	hsa-mir-151a	
QM073-2	F02	HmiRQP0125	MIMAT0000424	hsa-mir-128-1	
QM073-2	F03	HmiRQP0696	MIMAT0003260	hsa-mir-592	
QM073-2	F04	HmiRQP0479	MIMAT0000736	hsa-mir-381	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM073-2	F05	HmiRQP0197	MIMAT0002809	hsa-mir-146b	
QM073-2	F06	HmiRQP0761	MIMAT0003321	hsa-mir-651	
QM073-2	F07	HmiRQP0339	MIMAT0000279	hsa-mir-222	
QM073-2	F08	HmiRQP0451	MIMAT0000710	hsa-mir-365a	
QM073-2	F09	HmiRQP0617	MIMAT0003163	hsa-mir-539	
QM073-2	F10	HmiRQP0362	MIMAT0000085	hsa-mir-28	
QM073-2	F11	HmiRQP0786	MIMAT0000252	hsa-mir-7-1	
QM073-2	F12	HmiRQP0736	MIMAT0004809	hsa-mir-628	
QM073-2	G01	HmiRQP0242	MIMAT0006767	hsa-mir-1827	
QM073-2	G02	HmiRQP0779	MIMAT0003881	hsa-mir-668	
QM073-2	G03	HmiRQP0401	MIMAT0000089	hsa-mir-31	
QM073-2	G04	HmiRQP0278	MIMAT0002819	hsa-mir-193b	
QM073-2	G05	HmiRQP0156	MIMAT0000425	hsa-mir-130a	
QM073-2	G06	HmiRQP0854	MIMAT0000097	hsa-mir-99a	
QM073-2	G07	HmiRQP0509	MIMAT0001631	hsa-mir-451a	
QM073-2	G08	HmiRQP0018	MIMAT0000098	hsa-mir-100	
QM073-2	G09	HmiRQP0693	MIMAT0003258	hsa-mir-590	
QM073-2	G10	HmiRQP0832	MIMAT0000092	hsa-mir-92a-1	
QM073-2	G11	HmiRQP0421	MIMAT0000765	hsa-mir-335	
QM073-2	G12	HmiRQP0308	MIMAT0000462	hsa-mir-206	
QM073-2	H01	NC	NC	NC	
QM073-2	H02	NC	NC	NC	
QM073-2	H03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM073-2	H04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM073-2	H05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM073-2	H06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM073-2	H07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM073-2	H08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM073-2	H09	RT	RT	RT	
QM073-2	H10	RT	RT	RT	
QM073-2	H11	PCR	PCR	PCR	
QM073-2	H12	PCR	PCR	PCR	
QM073-3	A01	HmiRQP0276	MIMAT0000459	hsa-mir-193a	
QM073-3	A02	HmiRQP0263	MIMAT0000440	hsa-mir-191	
QM073-3	A03	HmiRQP0252	MIMAT0000457	hsa-mir-188	
QM073-3	A04	HmiRQP0248	MIMAT0000456	hsa-mir-186	
QM073-3	A05	HmiRQP0185	MIMAT0000433	hsa-mir-142	
QM073-3	A06	HmiRQP0181	MIMAT0000431	hsa-mir-140	
QM073-3	A07	HmiRQP0177	MIMAT0000430	hsa-mir-138-1	
QM073-3	A08	HmiRQP0168	MIMAT0000447	hsa-mir-134	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM073-3	A09	HmiRQP0491	MIMAT0001340	hsa-mir-423	
QM073-3	A10	HmiRQP0441	MIMAT0004676	hsa-mir-34b	
QM073-3	A11	HmiRQP0349	MIMAT0000080	hsa-mir-24-1	
QM073-3	A12	HmiRQP0319	MIMAT0000269	hsa-mir-212	
QM073-3	B01	HmiRQP0317	MIMAT0000267	hsa-mir-210	
QM073-3	B02	HmiRQP0309	MIMAT0000241	hsa-mir-208a	
QM073-3	B03	HmiRQP0227	MIMAT0000069	hsa-mir-16-1	
QM073-3	B04	HmiRQP0137	MIMAT0000242	hsa-mir-129-1	
QM073-3	B05	HmiRQP0373	MIMAT0000100	hsa-mir-29b-1	
QM073-3	B06	HmiRQP0354	MIMAT0000082	hsa-mir-26a-1	
QM073-3	B07	HmiRQP0332	MIMAT0000077	hsa-mir-22	
QM073-3	B08	HmiRQP0258	MIMAT0000458	hsa-mir-190a	
QM073-3	B09	HmiRQP0214	MIMAT0000439	hsa-mir-153-1	
QM073-3	B10	HmiRQP0175	MIMAT0000429	hsa-mir-137	
QM073-3	B11	HmiRQP0096	MIMAT0000423	hsa-mir-125b-1	
QM073-3	B12	HmiRQP0047	MIMAT0005865	hsa-mir-1202	
QM073-3	C01	HmiRQP0042	MIMAT0005798	hsa-mir-1185-1	
QM073-3	C02	HmiRQP0021	MIMAT0000099	hsa-mir-101-1	
QM073-3	C03	HmiRQP0044	MIMAT0000416	hsa-mir-1-1	
QM073-3	C04	HmiRQP0482	MIMAT0001075	hsa-mir-384	
QM073-3	C05	HmiRQP3590	MIMAT0027394	hsa-mir-6747	
QM073-3	C06	HmiRQP3602	MIMAT0027406	hsa-mir-6753	
QM073-3	C07	HmiRQP3604	MIMAT0027408	hsa-mir-6754	
QM073-3	C08	HmiRQP3620	MIMAT0027424	hsa-mir-6762	
QM073-3	C09	HmiRQP3670	MIMAT0027474	hsa-mir-6787	
QM073-3	C10	HmiRQP3778	MIMAT0027582	hsa-mir-6840	
QM073-3	C11	HmiRQP1317	MIMAT0012734	hsa-mir-711	
QM073-3	C12	HmiRQP0853	MIMAT0000096	hsa-mir-98	
QM073-3	D01	HmiRQP3618	MIMAT0027422	hsa-mir-6761	
QM073-3	E01	NC	NC	NC	
QM073-3	E02	NC	NC	NC	
QM073-3	E03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM073-3	E04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM073-3	E05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM073-3	E06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM073-3	E07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM073-3	E08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM073-3	E09	RT	RT	RT	
QM073-3	E10	RT	RT	RT	
QM073-3	E11	PCR	PCR	PCR	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM073-3	E12	PCR	PCR	PCR	