

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-01	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1	
QM004-01	A02	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM004-01	A03	HmiRQP0025	MIMAT0000102	hsa-miR-105-5p	
QM004-01	A04	HmiRQP0029	MIMAT0000680	hsa-miR-106b-5p	
QM004-01	A05	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM004-01	A06	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM004-01	A07	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM004-01	A08	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM004-01	A09	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM004-01	A10	HmiRQP0175	MIMAT0000429	hsa-miR-137	
QM004-01	A11	HmiRQP0180	MIMAT0000250	hsa-miR-139-5p	
QM004-01	A12	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM004-01	B01	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM004-01	B02	HmiRQP0214	MIMAT0000439	hsa-miR-153	
QM004-01	B03	HmiRQP0219	MIMAT0000452	hsa-miR-154-5p	
QM004-01	B04	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM004-01	B05	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM004-01	B06	HmiRQP0258	MIMAT0000458	hsa-miR-190a	
QM004-01	B07	HmiRQP0330	MIMAT0000276	hsa-miR-219-5p	
QM004-01	B08	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM004-01	B09	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM004-01	B10	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM004-01	B11	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM004-01	B12	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM004-01	C01	HmiRQP0857	MIMAT0000689	hsa-miR-99b-5p	
QM004-01	C02	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM004-01	C03	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM004-01	C04	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM004-01	C05	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM004-01	C06	HmiRQP0317	MIMAT0000267	hsa-miR-210	
QM004-01	C07	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM004-01	C08	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM004-01	C09	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM004-01	C10	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM004-01	C11	HmiRQP0459	MIMAT0000724	hsa-miR-372	
QM004-01	C12	HmiRQP0491	MIMAT0001340	hsa-miR-423-3p	
QM004-01	D01	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM004-01	D02	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM004-01	D03	HmiRQP0168	MIMAT0000447	hsa-miR-134	
QM004-01	D04	HmiRQP0173	MIMAT0000448	hsa-miR-136-5p	
QM004-01	D05	HmiRQP0177	MIMAT0000430	hsa-miR-138-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-01	D06	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM004-01	D07	HmiRQP0186	MIMAT0000434	hsa-miR-142-3p	
QM004-01	D08	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM004-01	D09	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM004-01	D10	HmiRQP0213	MIMAT0000438	hsa-miR-152	
QM004-01	D11	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM004-01	D12	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM004-01	E01	HmiRQP0251	MIMAT0000262	hsa-miR-187-3p	
QM004-01	E02	HmiRQP0252	MIMAT0000457	hsa-miR-188-5p	
QM004-01	E03	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM004-01	E04	HmiRQP0276	MIMAT0000459	hsa-miR-193a-3p	
QM004-01	E05	HmiRQP0280	MIMAT0000460	hsa-miR-194-5p	
QM004-01	E06	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM004-01	E07	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM004-01	E08	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM004-01	E09	HmiRQP0323	MIMAT0000273	hsa-miR-216a	
QM004-01	E10	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM004-01	E11	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM004-01	E12	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM004-01	F01	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM004-01	F02	HmiRQP0364	MIMAT0000690	hsa-miR-296-5p	
QM004-01	F03	HmiRQP0381	MIMAT0000684	hsa-miR-302a-3p	
QM004-01	F04	HmiRQP0404	MIMAT0000090	hsa-miR-32-5p	
QM004-01	F05	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM004-01	F06	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM004-01	F07	HmiRQP0461	MIMAT0000726	hsa-miR-373-3p	
QM004-01	F08	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM004-01	F09	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM004-01	F10	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM004-01	F11	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM004-01	F12	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM004-01	G01	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM004-01	G02	HmiRQP0325	MIMAT0000274	hsa-miR-217	
QM004-01	G03	HmiRQP0510	MIMAT0001635	hsa-miR-452-5p	
QM004-01	G04	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM004-01	G05	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM004-01	G06	HmiRQP0556	MIMAT0004778	hsa-miR-508-5p	
QM004-01	G07	HmiRQP0671	MIMAT0003237	hsa-miR-572	
QM004-01	G08	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM004-01	G09	HmiRQP0531	MIMAT0002807	hsa-miR-491-5p	
QM004-01	G10	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-01	G11	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM004-01	G12	HmiRQP0190	MIMAT0000436	hsa-miR-144-3p	
QM004-01	H01	NC	NC	NC	
QM004-01	H02	NC	NC	NC	
QM004-01	H03	HmiRQP9001	NR_002752	HK1	
QM004-01	H04	HmiRQP9011	NR_002750	HK2	
QM004-01	H05	HmiRQP9021	NR_002745	HK3	
QM004-01	H06	HmiRQP9051	NR_002746	HK4	
QM004-01	H07	HmiRQP9061	NR_002744	HK5	
QM004-01	H08	HmiRQP9071	NR_002450	HK6	
QM004-01	H09	RT	RT	RT	
QM004-01	H10	RT	RT	RT	
QM004-01	H11	PCR	PCR	PCR	
QM004-01	H12	PCR	PCR	PCR	
QM004-02	A01	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	
QM004-02	A02	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM004-02	A03	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM004-02	A04	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM004-02	A05	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM004-02	A06	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM004-02	A07	HmiRQP0389	MIMAT0005932	hsa-miR-302f	
QM004-02	A08	HmiRQP0540	MIMAT0002820	hsa-miR-497-5p	
QM004-02	A09	HmiRQP0569	MIMAT0002883	hsa-miR-514a-3p	
QM004-02	A10	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM004-02	A11	HmiRQP0388	MIMAT0005931	hsa-miR-302e	
QM004-02	A12	HmiRQP0485	MIMAT0002171	hsa-miR-410	
QM004-02	B01	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM004-02	B02	HmiRQP0262	MIMAT0004929	hsa-miR-190b	
QM004-02	B03	HmiRQP0486	MIMAT0003329	hsa-miR-411-5p	
QM004-02	B04	HmiRQP0736	MIMAT0004809	hsa-miR-628-5p	
QM004-02	B05	HmiRQP0661	MIMAT0003225	hsa-miR-561-3p	
QM004-02	B06	HmiRQP0434	MIMAT0004692	hsa-miR-340-5p	
QM004-02	B07	HmiRQP0735	MIMAT0003296	hsa-miR-627	
QM004-02	B08	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM004-02	B09	HmiRQP0476	MIMAT0000733	hsa-miR-379-5p	
QM004-02	B10	HmiRQP0596	MIMAT0002833	hsa-miR-520a-5p	
QM004-02	B11	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM004-02	B12	HmiRQP0413	MIMAT0000771	hsa-miR-325	
QM004-02	C01	HmiRQP0663	MIMAT0003227	hsa-miR-563	
QM004-02	C02	HmiRQP0334	MIMAT0000277	hsa-miR-220a	
QM004-02	C03	HmiRQP0566	MIMAT0002877	hsa-miR-513a-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-02	C04	HmiRQP0305	MIMAT0000264	hsa-miR-203	
QM004-02	C05	HmiRQP0362	MIMAT0000085	hsa-miR-28-5p	
QM004-02	C06	HmiRQP0416	MIMAT0001629	hsa-miR-329	
QM004-02	C07	HmiRQP0451	MIMAT0000710	hsa-miR-365a-3p	
QM004-02	C08	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM004-02	C09	HmiRQP0532	MIMAT0004765	hsa-miR-491-3p	
QM004-02	C10	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM004-02	C11	HmiRQP0212	MIMAT0004697	hsa-miR-151a-5p	
QM004-02	C12	HmiRQP0684	MIMAT0003247	hsa-miR-582-5p	
QM004-02	D01	HmiRQP0424	MIMAT0004695	hsa-miR-337-5p	
QM004-02	D02	HmiRQP0737	MIMAT0003297	hsa-miR-628-3p	
QM004-02	D03	HmiRQP0479	MIMAT0000736	hsa-miR-381	
QM004-02	D04	HmiRQP0574	MIMAT0006778	hsa-miR-516a-3p	hsa-miR-516b-3p
QM004-02	D05	HmiRQP0423	MIMAT0000754	hsa-miR-337-3p	
QM004-02	D06	HmiRQP0472	MIMAT0000730	hsa-miR-377-3p	
QM004-02	D07	HmiRQP0537	MIMAT0002817	hsa-miR-495	
QM004-02	D08	HmiRQP0125	MIMAT0000424	hsa-miR-128	
QM004-02	D09	HmiRQP0565	MIMAT0004777	hsa-miR-513a-3p	
QM004-02	D10	HmiRQP0683	MIMAT0004797	hsa-miR-582-3p	
QM004-02	D11	HmiRQP0211	MIMAT0000757	hsa-miR-151a-3p	
QM004-02	D12	HmiRQP0366	MIMAT0004450	hsa-miR-297	
QM004-02	E01	HmiRQP0410	MIMAT0000755	hsa-miR-323a-3p	
QM004-02	E02	HmiRQP0473	MIMAT0000732	hsa-miR-378a-3p	
QM004-02	E03	HmiRQP0369	MIMAT0002890	hsa-miR-299-5p	
QM004-02	E04	HmiRQP0562	MIMAT0002808	hsa-miR-511	
QM004-02	E05	HmiRQP0605	MIMAT0002854	hsa-miR-521	
QM004-02	E06	HmiRQP0446	MIMAT0000703	hsa-miR-361-5p	
QM004-02	E07	HmiRQP0481	MIMAT0000738	hsa-miR-383	
QM004-02	E08	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM004-02	E09	HmiRQP0557	MIMAT0002880	hsa-miR-508-3p	
QM004-02	E10	HmiRQP0610	MIMAT0002838	hsa-miR-525-5p	
QM004-02	E11	HmiRQP0758	MIMAT0003318	hsa-miR-648	
QM004-02	E12	HmiRQP0363	MIMAT0004502	hsa-miR-28-3p	
QM004-02	F01	HmiRQP0435	MIMAT0004694	hsa-miR-342-5p	
QM004-02	F02	HmiRQP0331	MIMAT0004675	hsa-miR-219-2-3p	
QM004-02	F03	HmiRQP0336	MIMAT0004915	hsa-miR-220c	
QM004-02	F04	HmiRQP0199	MIMAT0000251	hsa-miR-147a	
QM004-02	F05	HmiRQP0458	MIMAT0004687	hsa-miR-371a-5p	
QM004-02	F06	HmiRQP0452	MIMAT0000719	hsa-miR-367-3p	
QM004-02	F07	HmiRQP0850	MIMAT0000094	hsa-miR-95	
QM004-02	F08	HmiRQP0561	MIMAT0002882	hsa-miR-510	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-02	F09	HmiRQP0430	MIMAT0000091	hsa-miR-33a-5p	
QM004-02	F10	HmiRQP0516	MIMAT0003150	hsa-miR-455-5p	
QM004-02	F11	HmiRQP0712	MIMAT0003276	hsa-miR-608	
QM004-02	F12	HmiRQP0518	MIMAT0002173	hsa-miR-483-3p	
QM004-02	G01	HmiRQP0335	MIMAT0004908	hsa-miR-220b	
QM004-02	G02	HmiRQP0752	MIMAT0003312	hsa-miR-642a-5p	
QM004-02	G03	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM004-02	G04	HmiRQP0425	MIMAT0004701	hsa-miR-338-5p	
QM004-02	G05	HmiRQP0515	MIMAT0004784	hsa-miR-455-3p	
QM004-02	G06	HmiRQP0717	MIMAT0003281	hsa-miR-613	
QM004-02	G07	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM004-02	G08	HmiRQP0699	MIMAT0003263	hsa-miR-595	
QM004-02	G09	HmiRQP0528	MIMAT0002805	hsa-miR-489	
QM004-02	G10	HmiRQP0783	MIMAT0004926	hsa-miR-708-5p	
QM004-02	G11	HmiRQP0303	MIMAT0002811	hsa-miR-202-3p	
QM004-02	G12	HmiRQP0368	MIMAT0000687	hsa-miR-299-3p	
QM004-02	H01	NC	NC	NC	
QM004-02	H02	NC	NC	NC	
QM004-02	H03	HmiRQP9001	NR_002752	HK1	
QM004-02	H04	HmiRQP9011	NR_002750	HK2	
QM004-02	H05	HmiRQP9021	NR_002745	HK3	
QM004-02	H06	HmiRQP9051	NR_002746	HK4	
QM004-02	H07	HmiRQP9061	NR_002744	HK5	
QM004-02	H08	HmiRQP9071	NR_002450	HK6	
QM004-02	H09	RT	RT	RT	
QM004-02	H10	RT	RT	RT	
QM004-02	H11	PCR	PCR	PCR	
QM004-02	H12	PCR	PCR	PCR	
QM004-03	A01	HmiRQP0324	MIMAT0004959	hsa-miR-216b	
QM004-03	A02	HmiRQP0555	MIMAT0002879	hsa-miR-507	
QM004-03	A03	HmiRQP0615	MIMAT0002888	hsa-miR-532-5p	
QM004-03	A04	HmiRQP0490	MIMAT0001339	hsa-miR-422a	
QM004-03	A05	HmiRQP0536	MIMAT0002816	hsa-miR-494	
QM004-03	A06	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM004-03	A07	HmiRQP0534	MIMAT0003161	hsa-miR-493-3p	
QM004-03	A08	HmiRQP0728	MIMAT0003291	hsa-miR-622	
QM004-03	A09	HmiRQP0545	MIMAT0004773	hsa-miR-500a-5p	
QM004-03	A10	HmiRQP0572	MIMAT0004770	hsa-miR-516a-5p	
QM004-03	A11	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM004-03	A12	HmiRQP0237	MIMAT0002821	hsa-miR-181d	
QM004-03	B01	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-03	B02	HmiRQP0139	MIMAT0004605	hsa-miR-129-2-3p	
QM004-03	B03	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM004-03	B04	HmiRQP0412	MIMAT0000761	hsa-miR-324-5p	
QM004-03	B05	HmiRQP0686	MIMAT0003249	hsa-miR-584-5p	
QM004-03	B06	HmiRQP0426	MIMAT0000763	hsa-miR-338-3p	
QM004-03	B07	HmiRQP0480	MIMAT0000737	hsa-miR-382-5p	
QM004-03	B08	HmiRQP0546	MIMAT0002872	hsa-miR-501-5p	
QM004-03	B09	HmiRQP0419	MIMAT0004700	hsa-miR-331-5p	
QM004-03	B10	HmiRQP0488	MIMAT0002170	hsa-miR-412	
QM004-03	B11	HmiRQP0709	MIMAT0003273	hsa-miR-605	
QM004-03	B12	HmiRQP0570	MIMAT0002827	hsa-miR-515-3p	
QM004-03	C01	HmiRQP0581	MIMAT0002844	hsa-miR-518b	
QM004-03	C02	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM004-03	C03	HmiRQP0437	MIMAT0000772	hsa-miR-345-5p	
QM004-03	C04	HmiRQP0445	MIMAT0004682	hsa-miR-361-3p	
QM004-03	C05	HmiRQP0517	MIMAT0004761	hsa-miR-483-5p	
QM004-03	C06	HmiRQP0768	MIMAT0003335	hsa-miR-657	
QM004-03	C07	HmiRQP0457	MIMAT0000723	hsa-miR-371a-3p	
QM004-03	C08	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM004-03	C09	HmiRQP0495	MIMAT0003393	hsa-miR-425-5p	
QM004-03	C10	HmiRQP0522	MIMAT0004762	hsa-miR-486-3p	
QM004-03	C11	HmiRQP0533	MIMAT0002812	hsa-miR-492	
QM004-03	C12	HmiRQP0420	MIMAT0000760	hsa-miR-331-3p	
QM004-03	D01	HmiRQP0418	MIMAT0004693	hsa-miR-330-5p	
QM004-03	D02	HmiRQP0501	MIMAT0002814	hsa-miR-432-5p	
QM004-03	D03	HmiRQP0649	MIMAT0003214	hsa-miR-551a	
QM004-03	D04	HmiRQP0329	MIMAT0004567	hsa-miR-219-1-3p	
QM004-03	D05	HmiRQP0762	MIMAT0003322	hsa-miR-652-3p	
QM004-03	D06	HmiRQP0198	MIMAT0004766	hsa-miR-146b-3p	
QM004-03	D07	HmiRQP0436	MIMAT0000753	hsa-miR-342-3p	
QM004-03	D08	HmiRQP0492	MIMAT0004748	hsa-miR-423-5p	
QM004-03	D09	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM004-03	D10	HmiRQP0745	MIMAT0003305	hsa-miR-635	
QM004-03	D11	HmiRQP0811	MIMAT0004948	hsa-miR-885-3p	
QM004-03	D12	HmiRQP0520	MIMAT0002175	hsa-miR-485-5p	
QM004-03	E01	HmiRQP0414	MIMAT0000756	hsa-miR-326	
QM004-03	E02	HmiRQP0253	MIMAT0004613	hsa-miR-188-3p	
QM004-03	E03	HmiRQP0564	MIMAT0002822	hsa-miR-512-5p	
QM004-03	E04	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	
QM004-03	E05	HmiRQP0523	MIMAT0002177	hsa-miR-486-5p	
QM004-03	E06	HmiRQP0550	MIMAT0002874	hsa-miR-503	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-03	E07	HmiRQP0611	MIMAT0002839	hsa-miR-525-3p	
QM004-03	E08	HmiRQP0620	MIMAT0003340	hsa-miR-542-5p	
QM004-03	E09	HmiRQP0720	MIMAT0003283	hsa-miR-615-3p	
QM004-03	E10	HmiRQP0411	MIMAT0000762	hsa-miR-324-3p	
QM004-03	E11	HmiRQP0287	MIMAT0000227	hsa-miR-197-3p	
QM004-03	E12	HmiRQP0365	MIMAT0004679	hsa-miR-296-3p	
QM004-03	F01	HmiRQP0415	MIMAT0000752	hsa-miR-328	
QM004-03	F02	HmiRQP0179	MIMAT0004552	hsa-miR-139-3p	
QM004-03	F03	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM004-03	F04	HmiRQP0456	MIMAT0000722	hsa-miR-370	
QM004-03	F05	HmiRQP0718	MIMAT0003282	hsa-miR-614	
QM004-03	F06	HmiRQP0466	MIMAT0000728	hsa-miR-375	
QM004-03	F07	HmiRQP0417	MIMAT0000751	hsa-miR-330-3p	
QM004-03	F08	HmiRQP0277	MIMAT0004614	hsa-miR-193a-5p	
QM004-03	F09	HmiRQP0541	MIMAT0002824	hsa-miR-498	
QM004-03	F10	HmiRQP0438	MIMAT0000773	hsa-miR-346	
QM004-03	F11	HmiRQP0774	MIMAT0003326	hsa-miR-663a	
QM004-03	F12	HmiRQP0715	MIMAT0003279	hsa-miR-611	
QM004-03	G01	HmiRQP0747	MIMAT0003307	hsa-miR-637	
QM004-03	G02	HmiRQP0428	MIMAT0004702	hsa-miR-339-3p	
QM004-03	G03	HmiRQP0409	MIMAT0004696	hsa-miR-323a-5p	
QM004-03	G04	HmiRQP0772	MIMAT0003324	hsa-miR-661	
QM004-03	G05	HmiRQP0525	MIMAT0003180	hsa-miR-487b	
QM004-03	G06	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM004-03	G07	HmiRQP0463	MIMAT0000727	hsa-miR-374a-5p	
QM004-03	G08	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM004-03	G09	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM004-03	G10	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM004-03	G11	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	
QM004-03	G12	HmiRQP0204	MIMAT0000243	hsa-miR-148a-3p	
QM004-03	H01	NC	NC	NC	
QM004-03	H02	NC	NC	NC	
QM004-03	H03	HmiRQP9001	NR_002752	HK1	
QM004-03	H04	HmiRQP9011	NR_002750	HK2	
QM004-03	H05	HmiRQP9021	NR_002745	HK3	
QM004-03	H06	HmiRQP9051	NR_002746	HK4	
QM004-03	H07	HmiRQP9061	NR_002744	HK5	
QM004-03	H08	HmiRQP9071	NR_002450	HK6	
QM004-03	H09	RT	RT	RT	
QM004-03	H10	RT	RT	RT	
QM004-03	H11	PCR	PCR	PCR	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-03	H12	PCR	PCR	PCR	
QM004-04	A01	HmiRQP0469	MIMAT0002172	hsa-miR-376b	
QM004-04	A02	HmiRQP0467	MIMAT0000729	hsa-miR-376a-3p	
QM004-04	A03	HmiRQP0588	MIMAT0002869	hsa-miR-519a-3p	
QM004-04	A04	HmiRQP0470	MIMAT0000720	hsa-miR-376c	
QM004-04	A05	HmiRQP0853	MIMAT0000096	hsa-miR-98	
QM004-04	A06	HmiRQP0383	MIMAT0000715	hsa-miR-302b-3p	
QM004-04	A07	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM004-04	A08	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM004-04	A09	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM004-04	A10	HmiRQP0206	MIMAT0000759	hsa-miR-148b-3p	
QM004-04	A11	HmiRQP0322	MIMAT0000272	hsa-miR-215	
QM004-04	A12	HmiRQP0408	MIMAT0006764	hsa-miR-320d	
QM004-04	B01	HmiRQP0593	MIMAT0002829	hsa-miR-519e-3p	
QM004-04	B02	HmiRQP0387	MIMAT0000718	hsa-miR-302d-3p	
QM004-04	B03	HmiRQP0504	MIMAT0001541	hsa-miR-449a	
QM004-04	B04	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM004-04	B05	HmiRQP0314	MIMAT0001413	hsa-miR-20b-5p	
QM004-04	B06	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM004-04	B07	HmiRQP0378	MIMAT0000688	hsa-miR-301a-3p	
QM004-04	B08	HmiRQP0407	MIMAT0005793	hsa-miR-320c	
QM004-04	B09	HmiRQP0595	MIMAT0002834	hsa-miR-520a-3p	
QM004-04	B10	HmiRQP0592	MIMAT0002853	hsa-miR-519d	
QM004-04	B11	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM004-04	B12	HmiRQP0505	MIMAT0003327	hsa-miR-449b-5p	
QM004-04	C01	HmiRQP0379	MIMAT0004958	hsa-miR-301b	
QM004-04	C02	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM004-04	C03	HmiRQP0159	MIMAT0000691	hsa-miR-130b-3p	
QM004-04	C04	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM004-04	C05	HmiRQP0384	MIMAT0000717	hsa-miR-302c-3p	
QM004-04	C06	HmiRQP0318	MIMAT0000268	hsa-miR-211-5p	
QM004-04	C07	HmiRQP0547	MIMAT0004774	hsa-miR-501-3p	
QM004-04	C08	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM004-04	C09	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM004-04	C10	HmiRQP0006	MIMAT0000064	hsa-let-7c	
QM004-04	C11	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM004-04	C12	HmiRQP0170	MIMAT0000428	hsa-miR-135a-5p	
QM004-04	D01	HmiRQP0255	MIMAT0000072	hsa-miR-18a-5p	
QM004-04	D02	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM004-04	D03	HmiRQP0256	MIMAT0001412	hsa-miR-18b-5p	
QM004-04	D04	HmiRQP0598	MIMAT0002846	hsa-miR-520c-3p	



Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-04	D05	HmiRQP0172	MIMAT0000758	hsa-miR-135b-5p	
QM004-04	D06	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM004-04	D07	HmiRQP0575	MIMAT0002852	hsa-miR-517a-3p	
QM004-04	D08	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM004-04	D09	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM004-04	D10	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM004-04	D11	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM004-04	D12	HmiRQP0286	MIMAT0001080	hsa-miR-196b-5p	
QM004-04	E01	HmiRQP0601	MIMAT0002825	hsa-miR-520e	
QM004-04	E02	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM004-04	E03	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM004-04	E04	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM004-04	E05	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM004-04	E06	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM004-04	E07	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM004-04	E08	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM004-04	E09	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM004-04	E10	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM004-04	E11	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM004-04	E12	HmiRQP0405	MIMAT0000510	hsa-miR-320a	
QM004-04	F01	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM004-04	F02	HmiRQP0166	MIMAT0000427	hsa-miR-133a	
QM004-04	F03	HmiRQP1323	MIMAT0010251	hsa-miR-449c-5p	
QM004-04	F04	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	hsa-miR-199b-3p
QM004-04	F05	HmiRQP0612	MIMAT0005455	hsa-miR-520c-5p	hsa-miR-526a hsa-miR-518d-5p
QM004-04	F06	HmiRQP0462	MIMAT0004688	hsa-miR-374a-3p	
QM004-04	F07	HmiRQP0236	MIMAT0004559	hsa-miR-181c-3p	
QM004-04	F08	HmiRQP0218	MIMAT0000453	hsa-miR-154-3p	
QM004-04	F09	HmiRQP0471	MIMAT0004689	hsa-miR-377-5p	
QM004-04	F10	HmiRQP0500	MIMAT0002815	hsa-miR-432-3p	
QM004-04	F11	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM004-04	F12	HmiRQP0275	MIMAT0004543	hsa-miR-192-3p	
QM004-04	G01	HmiRQP0233	MIMAT0004558	hsa-miR-181a-2-3p	
QM004-04	G02	HmiRQP0176	MIMAT0004607	hsa-miR-138-1-3p	
QM004-04	G03	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM004-04	G04	HmiRQP0171	MIMAT0004698	hsa-miR-135b-3p	
QM004-04	G05	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM004-04	G06	HmiRQP0299	MIMAT0004571	hsa-miR-200b-5p	
QM004-04	G07	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM004-04	G08	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-04	G09	HmiRQP0005	MIMAT0004483	hsa-let-7c*	
QM004-04	G10	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM004-04	G11	HmiRQP0009	MIMAT0004485	hsa-let-7e-3p	
QM004-04	G12	HmiRQP0011	MIMAT0004486	hsa-let-7f-1-3p	
QM004-04	H01	NC	NC	NC	
QM004-04	H02	NC	NC	NC	
QM004-04	H03	HmiRQP9001	NR_002752	HK1	
QM004-04	H04	HmiRQP9011	NR_002750	HK2	
QM004-04	H05	HmiRQP9021	NR_002745	HK3	
QM004-04	H06	HmiRQP9051	NR_002746	HK4	
QM004-04	H07	HmiRQP9061	NR_002744	HK5	
QM004-04	H08	HmiRQP9071	NR_002450	HK6	
QM004-04	H09	RT	RT	RT	
QM004-04	H10	RT	RT	RT	
QM004-04	H11	PCR	PCR	PCR	
QM004-04	H12	PCR	PCR	PCR	
QM004-05	A01	HmiRQP0013	MIMAT0004487	hsa-let-7f-2-3p	
QM004-05	A02	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM004-05	A03	HmiRQP0019	MIMAT0004512	hsa-miR-100-3p	
QM004-05	A04	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM004-05	A05	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM004-05	A06	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM004-05	A07	HmiRQP0033	MIMAT0004556	hsa-miR-10b-3p	
QM004-05	A08	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM004-05	A09	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM004-05	A10	HmiRQP0136	MIMAT0004548	hsa-miR-129-1-3p	
QM004-05	A11	HmiRQP0158	MIMAT0004680	hsa-miR-130b-5p	
QM004-05	A12	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM004-05	B01	HmiRQP0169	MIMAT0004595	hsa-miR-135a-3p	
QM004-05	B02	HmiRQP0178	MIMAT0004596	hsa-miR-138-2-3p	
QM004-05	B03	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM004-05	B04	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM004-05	B05	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM004-05	B06	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM004-05	B07	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM004-05	B08	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM004-05	B09	HmiRQP0249	MIMAT0004612	hsa-miR-186-3p	
QM004-05	B10	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM004-05	B11	HmiRQP0304	MIMAT0002810	hsa-miR-202-5p	
QM004-05	B12	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM004-05	C01	HmiRQP0321	MIMAT0004564	hsa-miR-214-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-05	C02	HmiRQP0328	MIMAT0004566	hsa-miR-218-2-3p	
QM004-05	C03	HmiRQP0333	MIMAT0004495	hsa-miR-22-5p	
QM004-05	C04	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM004-05	C05	HmiRQP0345	MIMAT0004496	hsa-miR-23a-5p	
QM004-05	C06	HmiRQP0347	MIMAT0004587	hsa-miR-23b-5p	
QM004-05	C07	HmiRQP0348	MIMAT0000079	hsa-miR-24-1-5p	
QM004-05	C08	HmiRQP0350	MIMAT0004497	hsa-miR-24-2-5p	
QM004-05	C09	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM004-05	C10	HmiRQP0353	MIMAT0004499	hsa-miR-26a-1-3p	
QM004-05	C11	HmiRQP0355	MIMAT0004681	hsa-miR-26a-2-3p	
QM004-05	C12	HmiRQP0360	MIMAT0004588	hsa-miR-27b-5p	
QM004-05	D01	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM004-05	D02	HmiRQP0372	MIMAT0004514	hsa-miR-29b-1-5p	
QM004-05	D03	HmiRQP0374	MIMAT0004515	hsa-miR-29b-2-5p	
QM004-05	D04	HmiRQP0376	MIMAT0004673	hsa-miR-29c-5p	
QM004-05	D05	HmiRQP0394	MIMAT0004674	hsa-miR-30c-1-3p	
QM004-05	D06	HmiRQP0397	MIMAT0004551	hsa-miR-30d-3p	
QM004-05	D07	HmiRQP0402	MIMAT0004504	hsa-miR-31-3p	
QM004-05	D08	HmiRQP0403	MIMAT0004505	hsa-miR-32-3p	
QM004-05	D09	HmiRQP0422	MIMAT0004703	hsa-miR-335-3p	
QM004-05	D10	HmiRQP0429	MIMAT0004506	hsa-miR-33a-3p	
QM004-05	D11	HmiRQP0433	MIMAT0000750	hsa-miR-340-3p	
QM004-05	D12	HmiRQP0442	MIMAT0000685	hsa-miR-34b-5p	
QM004-05	E01	HmiRQP0453	MIMAT0004686	hsa-miR-367-5p	
QM004-05	E02	HmiRQP0465	MIMAT0004956	hsa-miR-374b-3p	
QM004-05	E03	HmiRQP0474	MIMAT0000731	hsa-miR-378a-5p	
QM004-05	E04	HmiRQP0493	MIMAT0004749	hsa-miR-424-3p	
QM004-05	E05	HmiRQP0539	MIMAT0004768	hsa-miR-497-3p	
QM004-05	E06	HmiRQP0544	MIMAT0002871	hsa-miR-500a-3p	
QM004-05	E07	HmiRQP0576	MIMAT0002851	hsa-miR-517-5p	
QM004-05	E08	HmiRQP0784	MIMAT0004927	hsa-miR-708-3p	
QM004-05	E09	HmiRQP0785	MIMAT0004553	hsa-miR-7-1-3p	
QM004-05	E10	HmiRQP0787	MIMAT0004554	hsa-miR-7-2-3p	
QM004-05	E11	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM004-05	E12	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM004-05	F01	HmiRQP0855	MIMAT0004511	hsa-miR-99a-3p	
QM004-05	F02	HmiRQP0856	MIMAT0004678	hsa-miR-99b-3p	
QM004-05	F03	HmiRQP1295	MIMAT0009198	hsa-miR-224-3p	
QM004-05	F04	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM004-05	F05	HmiRQP1301	MIMAT0009203	hsa-miR-449b-3p	
QM004-05	F06	HmiRQP1309	MIMAT0010195	hsa-let-7a-2-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM004-05	F07	HmiRQP1523	MIMAT0015072	hsa-miR-320e	
QM004-05	F08	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM004-05	F09	HmiRQP0097	MIMAT0004603	hsa-miR-125b-2-3p	
QM004-05	F10	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM004-05	F11	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM004-05	F12	HmiRQP0205	MIMAT0004699	hsa-miR-148b-5p	
QM004-05	G01	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM004-05	G02	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM004-05	G03	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM004-05	G04	HmiRQP0224	MIMAT0004586	hsa-miR-15b-3p	
QM004-05	G05	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM004-05	G06	HmiRQP0228	MIMAT0004518	hsa-miR-16-2-3p	
QM004-05	G07	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM004-05	G08	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM004-05	G09	HmiRQP0243	MIMAT0004560	hsa-miR-183-3p	
QM004-05	G10	HmiRQP0257	MIMAT0004751	hsa-miR-18b-3p	
QM004-05	G11	HmiRQP0264	MIMAT0001618	hsa-miR-191-3p	
QM004-05	G12	HmiRQP0279	MIMAT0004767	hsa-miR-193b-5p	
QM004-05	H01	NC	NC	NC	
QM004-05	H02	NC	NC	NC	
QM004-05	H03	HmiRQP9001	NR_002752	HK1	
QM004-05	H04	HmiRQP9011	NR_002750	HK2	
QM004-05	H05	HmiRQP9021	NR_002745	HK3	
QM004-05	H06	HmiRQP9051	NR_002746	HK4	
QM004-05	H07	HmiRQP9061	NR_002744	HK5	
QM004-05	H08	HmiRQP9071	NR_002450	HK6	
QM004-05	H09	RT	RT	RT	
QM004-05	H10	RT	RT	RT	
QM004-05	H11	PCR	PCR	PCR	
QM004-05	H12	PCR	PCR	PCR	