

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM013-01	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1	
QM013-01	A02	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM013-01	A03	HmiRQP0025	MIMAT0000102	hsa-miR-105-5p	
QM013-01	A04	HmiRQP0029	MIMAT0000680	hsa-miR-106b-5p	
QM013-01	A05	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM013-01	A06	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM013-01	A07	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM013-01	A08	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM013-01	A09	HmiRQP0175	MIMAT0000429	hsa-miR-137	
QM013-01	A10	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM013-01	A11	HmiRQP0219	MIMAT0000452	hsa-miR-154-5p	
QM013-01	A12	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM013-01	B01	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM013-01	B02	HmiRQP0258	MIMAT0000458	hsa-miR-190a	
QM013-01	B03	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM013-01	B04	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM013-01	B05	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM013-01	B06	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM013-01	B07	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM013-01	B08	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM013-01	B09	HmiRQP0317	MIMAT0000267	hsa-miR-210	
QM013-01	B10	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM013-01	B11	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM013-01	B12	HmiRQP0168	MIMAT0000447	hsa-miR-134	
QM013-01	C01	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM013-01	C02	HmiRQP0186	MIMAT0000434	hsa-miR-142-3p	
QM013-01	C03	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM013-01	C04	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM013-01	C05	HmiRQP0213	MIMAT0000438	hsa-miR-152	
QM013-01	C06	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM013-01	C07	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM013-01	C08	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM013-01	C09	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM013-01	C10	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM013-01	C11	HmiRQP0364	MIMAT0000690	hsa-miR-296-5p	
QM013-01	C12	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM013-01	D01	HmiRQP0461	MIMAT0000726	hsa-miR-373-3p	
QM013-01	D02	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM013-01	D03	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM013-01	D04	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM013-01	D05	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM013-01	D06	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM013-01	D07	HmiRQP0556	MIMAT0004778	hsa-miR-508-5p	
QM013-01	D08	HmiRQP0671	MIMAT0003237	hsa-miR-572	
QM013-01	D09	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM013-01	D10	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM013-01	D11	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	
QM013-01	D12	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM013-01	E01	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM013-01	E02	HmiRQP0569	MIMAT0002883	hsa-miR-514a-3p	
QM013-01	E03	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM013-01	E04	HmiRQP0485	MIMAT0002171	hsa-miR-410	
QM013-01	E05	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM013-01	E06	HmiRQP0735	MIMAT0003296	hsa-miR-627	
QM013-01	E07	HmiRQP0476	MIMAT0000733	hsa-miR-379-5p	
QM013-01	E08	HmiRQP0596	MIMAT0002833	hsa-miR-520a-5p	
QM013-01	E09	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM013-01	E10	HmiRQP0413	MIMAT0000771	hsa-miR-325	
QM013-01	E11	HmiRQP0305	MIMAT0000264	hsa-miR-203	
QM013-01	E12	HmiRQP0451	MIMAT0000710	hsa-miR-365a-3p	
QM013-01	F01	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM013-01	F02	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM013-01	F03	HmiRQP0424	MIMAT0004695	hsa-miR-337-5p	
QM013-01	F04	HmiRQP0479	MIMAT0000736	hsa-miR-381	
QM013-01	F05	HmiRQP0423	MIMAT0000754	hsa-miR-337-3p	
QM013-01	F06	HmiRQP0472	MIMAT0000730	hsa-miR-377-3p	
QM013-01	F07	HmiRQP0537	MIMAT0002817	hsa-miR-495	
QM013-01	F08	HmiRQP0473	MIMAT0000732	hsa-miR-378a-3p	
QM013-01	F09	HmiRQP0446	MIMAT0000703	hsa-miR-361-5p	
QM013-01	F10	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM013-01	F11	HmiRQP0557	MIMAT0002880	hsa-miR-508-3p	
QM013-01	F12	HmiRQP0758	MIMAT0003318	hsa-miR-648	
QM013-01	G01	HmiRQP0199	MIMAT0000251	hsa-miR-147a	
QM013-01	G02	HmiRQP0712	MIMAT0003276	hsa-miR-608	
QM013-01	G03	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM013-01	G04	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM013-01	G05	HmiRQP0555	MIMAT0002879	hsa-miR-507	
QM013-01	G06	HmiRQP0490	MIMAT0001339	hsa-miR-422a	
QM013-01	G07	HmiRQP0536	MIMAT0002816	hsa-miR-494	
QM013-01	G08	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM013-01	G09	HmiRQP0534	MIMAT0003161	hsa-miR-493-3p	
QM013-01	G10	HmiRQP0728	MIMAT0003291	hsa-miR-622	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM013-01	G11	HmiRQP0572	MIMAT0004770	hsa-miR-516a-5p	
QM013-01	G12	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM013-01	H01	NC	NC	NC	
QM013-01	H02	NC	NC	NC	
QM013-01	H03	HmiRQP9001	NR_002752	HK1	
QM013-01	H04	HmiRQP9011	NR_002750	HK2	
QM013-01	H05	HmiRQP9021	NR_002745	HK3	
QM013-01	H06	HmiRQP9051	NR_002746	HK4	
QM013-01	H07	HmiRQP9061	NR_002744	HK5	
QM013-01	H08	HmiRQP9071	NR_002450	HK6	
QM013-01	H09	RT	RT	RT	
QM013-01	H10	RT	RT	RT	
QM013-01	H11	PCR	PCR	PCR	
QM013-01	H12	PCR	PCR	PCR	
QM013-02	A01	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM013-02	A02	HmiRQP0139	MIMAT0004605	hsa-miR-129-2-3p	
QM013-02	A03	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM013-02	A04	HmiRQP0419	MIMAT0004700	hsa-miR-331-5p	
QM013-02	A05	HmiRQP0709	MIMAT0003273	hsa-miR-605	
QM013-02	A06	HmiRQP0570	MIMAT0002827	hsa-miR-515-3p	
QM013-02	A07	HmiRQP0581	MIMAT0002844	hsa-miR-518b	
QM013-02	A08	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM013-02	A09	HmiRQP0445	MIMAT0004682	hsa-miR-361-3p	
QM013-02	A10	HmiRQP0517	MIMAT0004761	hsa-miR-483-5p	
QM013-02	A11	HmiRQP0768	MIMAT0003335	hsa-miR-657	
QM013-02	A12	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM013-02	B01	HmiRQP0533	MIMAT0002812	hsa-miR-492	
QM013-02	B02	HmiRQP0420	MIMAT0000760	hsa-miR-331-3p	
QM013-02	B03	HmiRQP0501	MIMAT0002814	hsa-miR-432-5p	
QM013-02	B04	HmiRQP0649	MIMAT0003214	hsa-miR-551a	
QM013-02	B05	HmiRQP0198	MIMAT0004766	hsa-miR-146b-3p	
QM013-02	B06	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM013-02	B07	HmiRQP0745	MIMAT0003305	hsa-miR-635	
QM013-02	B08	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	
QM013-02	B09	HmiRQP0365	MIMAT0004679	hsa-miR-296-3p	
QM013-02	B10	HmiRQP0456	MIMAT0000722	hsa-miR-370	
QM013-02	B11	HmiRQP0466	MIMAT0000728	hsa-miR-375	
QM013-02	B12	HmiRQP0541	MIMAT0002824	hsa-miR-498	
QM013-02	C01	HmiRQP0438	MIMAT0000773	hsa-miR-346	
QM013-02	C02	HmiRQP0774	MIMAT0003326	hsa-miR-663a	
QM013-02	C03	HmiRQP0715	MIMAT0003279	hsa-miR-611	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM013-02	C04	HmiRQP0747	MIMAT0003307	hsa-miR-637	
QM013-02	C05	HmiRQP0525	MIMAT0003180	hsa-miR-487b	
QM013-02	C06	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM013-02	C07	HmiRQP0463	MIMAT0000727	hsa-miR-374a-5p	
QM013-02	C08	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM013-02	C09	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM013-02	C10	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM013-02	C11	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	
QM013-02	C12	HmiRQP0469	MIMAT0002172	hsa-miR-376b	
QM013-02	D01	HmiRQP0467	MIMAT0000729	hsa-miR-376a-3p	
QM013-02	D02	HmiRQP0588	MIMAT0002869	hsa-miR-519a-3p	
QM013-02	D03	HmiRQP0853	MIMAT0000096	hsa-miR-98	
QM013-02	D04	HmiRQP0383	MIMAT0000715	hsa-miR-302b-3p	
QM013-02	D05	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM013-02	D06	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM013-02	D07	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM013-02	D08	HmiRQP0593	MIMAT0002829	hsa-miR-519e-3p	
QM013-02	D09	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM013-02	D10	HmiRQP0595	MIMAT0002834	hsa-miR-520a-3p	
QM013-02	D11	HmiRQP0592	MIMAT0002853	hsa-miR-519d	
QM013-02	D12	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM013-02	E01	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM013-02	E02	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM013-02	E03	HmiRQP0384	MIMAT0000717	hsa-miR-302c-3p	
QM013-02	E04	HmiRQP0318	MIMAT0000268	hsa-miR-211-5p	
QM013-02	E05	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM013-02	E06	HmiRQP0006	MIMAT0000064	hsa-let-7c	
QM013-02	E07	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM013-02	E08	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM013-02	E09	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM013-02	E10	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM013-02	E11	HmiRQP0601	MIMAT0002825	hsa-miR-520e	
QM013-02	E12	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM013-02	F01	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM013-02	F02	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM013-02	F03	HmiRQP0166	MIMAT0000427	hsa-miR-133a	
QM013-02	F04	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM013-02	F05	HmiRQP0462	MIMAT0004688	hsa-miR-374a-3p	
QM013-02	F06	HmiRQP0218	MIMAT0000453	hsa-miR-154-3p	
QM013-02	F07	HmiRQP0471	MIMAT0004689	hsa-miR-377-5p	
QM013-02	F08	HmiRQP0500	MIMAT0002815	hsa-miR-432-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM013-02	F09	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM013-02	F10	HmiRQP0299	MIMAT0004571	hsa-miR-200b-5p	
QM013-02	F11	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM013-02	F12	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM013-02	G01	HmiRQP0005	MIMAT0004483	hsa-let-7c*	
QM013-02	G02	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM013-02	G03	HmiRQP0009	MIMAT0004485	hsa-let-7e-3p	
QM013-02	G04	HmiRQP0011	MIMAT0004486	hsa-let-7f-1-3p	
QM013-02	G05	HmiRQP0013	MIMAT0004487	hsa-let-7f-2-3p	
QM013-02	G06	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM013-02	G07	HmiRQP0019	MIMAT0004512	hsa-miR-100-3p	
QM013-02	G08	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM013-02	G09	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM013-02	G10	HmiRQP0033	MIMAT0004556	hsa-miR-10b-3p	
QM013-02	G11	HmiRQP0136	MIMAT0004548	hsa-miR-129-1-3p	
QM013-02	G12	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM013-02	H01	NC	NC	NC	
QM013-02	H02	NC	NC	NC	
QM013-02	H03	HmiRQP9001	NR_002752	HK1	
QM013-02	H04	HmiRQP9011	NR_002750	HK2	
QM013-02	H05	HmiRQP9021	NR_002745	HK3	
QM013-02	H06	HmiRQP9051	NR_002746	HK4	
QM013-02	H07	HmiRQP9061	NR_002744	HK5	
QM013-02	H08	HmiRQP9071	NR_002450	HK6	
QM013-02	H09	RT	RT	RT	
QM013-02	H10	RT	RT	RT	
QM013-02	H11	PCR	PCR	PCR	
QM013-02	H12	PCR	PCR	PCR	