

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM037-01	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1	
QM037-01	A02	HmiRQP0029	MIMAT0000680	hsa-miR-106b-5p	
QM037-01	A03	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM037-01	A04	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM037-01	A05	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM037-01	A06	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM037-01	A07	HmiRQP0180	MIMAT0000250	hsa-miR-139-5p	
QM037-01	A08	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM037-01	A09	HmiRQP0219	MIMAT0000452	hsa-miR-154-5p	
QM037-01	A10	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM037-01	A11	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM037-01	A12	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM037-01	B01	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM037-01	B02	HmiRQP0857	MIMAT0000689	hsa-miR-99b-5p	
QM037-01	B03	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM037-01	B04	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM037-01	B05	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM037-01	B06	HmiRQP0459	MIMAT0000724	hsa-miR-372	
QM037-01	B07	HmiRQP0491	MIMAT0001340	hsa-miR-423-3p	
QM037-01	B08	HmiRQP0168	MIMAT0000447	hsa-miR-134	
QM037-01	B09	HmiRQP0173	MIMAT0000448	hsa-miR-136-5p	
QM037-01	B10	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM037-01	B11	HmiRQP0186	MIMAT0000434	hsa-miR-142-3p	
QM037-01	B12	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM037-01	C01	HmiRQP0213	MIMAT0000438	hsa-miR-152	
QM037-01	C02	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM037-01	C03	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM037-01	C04	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM037-01	C05	HmiRQP0280	MIMAT0000460	hsa-miR-194-5p	
QM037-01	C06	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM037-01	C07	HmiRQP0364	MIMAT0000690	hsa-miR-296-5p	
QM037-01	C08	HmiRQP0404	MIMAT0000090	hsa-miR-32-5p	
QM037-01	C09	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM037-01	C10	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM037-01	C11	HmiRQP0543	MIMAT0002870	hsa-miR-499a-5p	
QM037-01	C12	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM037-01	D01	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM037-01	D02	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM037-01	D03	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM037-01	D04	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM037-01	D05	HmiRQP0693	MIMAT0003258	hsa-miR-590-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM037-01	D06	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM037-01	D07	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM037-01	D08	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM037-01	D09	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM037-01	D10	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM037-01	D11	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM037-01	D12	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM037-01	E01	HmiRQP0540	MIMAT0002820	hsa-miR-497-5p	
QM037-01	E02	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM037-01	E03	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM037-01	E04	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM037-01	E05	HmiRQP0362	MIMAT0000085	hsa-miR-28-5p	
QM037-01	E06	HmiRQP0212	MIMAT0004697	hsa-miR-151a-5p	
QM037-01	E07	HmiRQP0537	MIMAT0002817	hsa-miR-495	
QM037-01	E08	HmiRQP0125	MIMAT0000424	hsa-miR-128	
QM037-01	E09	HmiRQP0211	MIMAT0000757	hsa-miR-151-3p	
QM037-01	E10	HmiRQP0366	MIMAT0004450	hsa-miR-297	
QM037-01	E11	HmiRQP0521	MIMAT0002176	hsa-miR-485-3p	
QM037-01	E12	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM037-01	F01	HmiRQP0812	MIMAT0004947	hsa-miR-885-5p	
QM037-01	F02	HmiRQP0363	MIMAT0004502	hsa-miR-28-3p	
QM037-01	F03	HmiRQP0739	MIMAT0004810	hsa-miR-629-5p	
QM037-01	F04	HmiRQP0850	MIMAT0000094	hsa-miR-95	
QM037-01	F05	HmiRQP0430	MIMAT0000091	hsa-miR-33a-5p	
QM037-01	F06	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM037-01	F07	HmiRQP0615	MIMAT0002888	hsa-miR-532-5p	
QM037-01	F08	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM037-01	F09	HmiRQP0771	MIMAT0003338	hsa-miR-660-5p	
QM037-01	F10	HmiRQP0545	MIMAT0004773	hsa-miR-500a-5p	
QM037-01	F11	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM037-01	F12	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM037-01	G01	HmiRQP0412	MIMAT0000761	hsa-miR-324-5p	
QM037-01	G02	HmiRQP0686	MIMAT0003249	hsa-miR-584-5p	
QM037-01	G03	HmiRQP0426	MIMAT0000763	hsa-miR-338-3p	
QM037-01	G04	HmiRQP0480	MIMAT0000737	hsa-miR-382-5p	
QM037-01	G05	HmiRQP0622	MIMAT0004954	hsa-miR-543	
QM037-01	G06	HmiRQP0709	MIMAT0003273	hsa-miR-605	
QM037-01	G07	HmiRQP0445	MIMAT0004682	hsa-miR-361-3p	
QM037-01	G08	HmiRQP0484	MIMAT0001639	hsa-miR-409-3p	
QM037-01	G09	HmiRQP0495	MIMAT0003393	hsa-miR-425-5p	
QM037-01	G10	HmiRQP0489	MIMAT0003339	hsa-miR-421	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM037-01	G11	HmiRQP0420	MIMAT0000760	hsa-miR-331-3p	
QM037-01	G12	HmiRQP0436	MIMAT0000753	hsa-miR-342-3p	
QM037-01	H01	NC	NC	NC	
QM037-01	H02	NC	NC	NC	
QM037-01	H03	HmiRQP9001	NR_002752	HK1	
QM037-01	H04	HmiRQP9011	NR_002750	HK2	
QM037-01	H05	HmiRQP9021	NR_002745	HK3	
QM037-01	H06	HmiRQP9051	NR_002746	HK4	
QM037-01	H07	HmiRQP9061	NR_002744	HK5	
QM037-01	H08	HmiRQP9071	NR_002450	HK6	
QM037-01	H09	RT	RT	RT	
QM037-01	H10	RT	RT	RT	
QM037-01	H11	PCR	PCR	PCR	
QM037-01	H12	PCR	PCR	PCR	
QM037-02	A01	HmiRQP0492	MIMAT0004748	hsa-miR-423-5p	
QM037-02	A02	HmiRQP0414	MIMAT0000756	hsa-miR-326	
QM037-02	A03	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	
QM037-02	A04	HmiRQP0523	MIMAT0002177	hsa-miR-486-5p	
QM037-02	A05	HmiRQP0519	MIMAT0002174	hsa-miR-484	
QM037-02	A06	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM037-02	A07	HmiRQP0411	MIMAT0000762	hsa-miR-324-3p	
QM037-02	A08	HmiRQP0415	MIMAT0000752	hsa-miR-328	
QM037-02	A09	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM037-02	A10	HmiRQP0466	MIMAT0000728	hsa-miR-375	
QM037-02	A11	HmiRQP0277	MIMAT0004614	hsa-miR-193a-5p	
QM037-02	A12	HmiRQP0438	MIMAT0000773	hsa-miR-346	
QM037-02	B01	HmiRQP0428	MIMAT0004702	hsa-miR-339-3p	
QM037-02	B02	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM037-02	B03	HmiRQP0463	MIMAT0000727	hsa-miR-374a-5p	
QM037-02	B04	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM037-02	B05	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM037-02	B06	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM037-02	B07	HmiRQP0470	MIMAT0000720	hsa-miR-376c	
QM037-02	B08	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM037-02	B09	HmiRQP0314	MIMAT0001413	hsa-miR-20b-5p	
QM037-02	B10	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM037-02	B11	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM037-02	B12	HmiRQP0318	MIMAT0000268	hsa-miR-211-5p	
QM037-02	C01	HmiRQP0548	MIMAT0004775	hsa-miR-502-3p	
QM037-02	C02	HmiRQP0547	MIMAT0004774	hsa-miR-501-3p	
QM037-02	C03	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM037-02	C04	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM037-02	C05	HmiRQP0006	MIMAT0000064	hsa-let-7c	
QM037-02	C06	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM037-02	C07	HmiRQP0256	MIMAT0001412	hsa-miR-18b-5p	
QM037-02	C08	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM037-02	C09	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM037-02	C10	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM037-02	C11	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM037-02	C12	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM037-02	D01	HmiRQP0405	MIMAT0000510	hsa-miR-320a	
QM037-02	D02	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM037-02	D03	HmiRQP0166	MIMAT0000427	hsa-miR-133a	
QM037-02	D04	HmiRQP1329	MIMAT0009449	hsa-miR-1974	
QM037-02	D05	HmiRQP1333	MIMAT0010133	hsa-miR-2110	
QM037-02	D06	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	hsa-miR-199b-3p
QM037-02	D07	HmiRQP0468	MIMAT0003386	hsa-miR-376a-5p	
QM037-02	D08	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM037-02	D09	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM037-02	D10	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM037-02	D11	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM037-02	D12	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM037-02	E01	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM037-02	E02	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM037-02	E03	HmiRQP0158	MIMAT0004680	hsa-miR-130b-5p	
QM037-02	E04	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM037-02	E05	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM037-02	E06	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM037-02	E07	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM037-02	E08	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM037-02	E09	HmiRQP0333	MIMAT0004495	hsa-miR-22-5p	
QM037-02	E10	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM037-02	E11	HmiRQP0345	MIMAT0004496	hsa-miR-23a-5p	
QM037-02	E12	HmiRQP0347	MIMAT0004587	hsa-miR-23b-5p	
QM037-02	F01	HmiRQP0348	MIMAT0000079	hsa-miR-24-1-5p	
QM037-02	F02	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM037-02	F03	HmiRQP0360	MIMAT0004588	hsa-miR-27b-5p	
QM037-02	F04	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM037-02	F05	HmiRQP0374	MIMAT0004515	hsa-miR-29b-2-5p	
QM037-02	F06	HmiRQP0376	MIMAT0004673	hsa-miR-29c-5p	
QM037-02	F07	HmiRQP0474	MIMAT0000731	hsa-miR-378a-5p	
QM037-02	F08	HmiRQP0553	MIMAT0004776	hsa-miR-505-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM037-02	F09	HmiRQP0650	MIMAT0004794	hsa-miR-551b-5p	
QM037-02	F10	HmiRQP0834	MIMAT0004792	hsa-miR-92b-5p	
QM037-02	F11	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM037-02	F12	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM037-02	G01	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM037-02	G02	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM037-02	G03	HmiRQP0205	MIMAT0004699	hsa-miR-148b-5p	
QM037-02	G04	HmiRQP0224	MIMAT0004586	hsa-miR-15b-3p	
QM037-02	G05	HmiRQP0228	MIMAT0004518	hsa-miR-16-2-3p	
QM037-02	G06	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM037-02	G07	HmiRQP0279	MIMAT0004767	hsa-miR-193b-5p	
QM037-02	G08	HmiRQP0340	MIMAT0004569	hsa-miR-222-5p	
QM037-02	G09	HmiRQP0358	MIMAT0004501	hsa-miR-27a-5p	
QM037-02	G10	HmiRQP0450	MIMAT0003385	hsa-miR-363-5p	
QM037-02	G11	HmiRQP0496	MIMAT0001343	hsa-miR-425-3p	
QM037-02	G12	HmiRQP0831	MIMAT0004507	hsa-miR-92a-1-5p	
QM037-02	H01	NC	NC	NC	
QM037-02	H02	NC	NC	NC	
QM037-02	H03	HmiRQP9001	NR_002752	HK1	
QM037-02	H04	HmiRQP9011	NR_002750	HK2	
QM037-02	H05	HmiRQP9021	NR_002745	HK3	
QM037-02	H06	HmiRQP9051	NR_002746	HK4	
QM037-02	H07	HmiRQP9061	NR_002744	HK5	
QM037-02	H08	HmiRQP9071	NR_002450	HK6	
QM037-02	H09	RT	RT	RT	
QM037-02	H10	RT	RT	RT	
QM037-02	H11	PCR	PCR	PCR	
QM037-02	H12	PCR	PCR	PCR	