

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM035-01	A01	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM035-01	A02	HmiRQP0219	MIMAT0000452	hsa-miR-154-5p	
QM035-01	A03	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM035-01	A04	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM035-01	A05	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM035-01	A06	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM035-01	A07	HmiRQP0857	MIMAT0000689	hsa-miR-99b-5p	
QM035-01	A08	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM035-01	A09	HmiRQP0317	MIMAT0000267	hsa-miR-210	
QM035-01	A10	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM035-01	A11	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM035-01	A12	HmiRQP0491	MIMAT0001340	hsa-miR-423-3p	
QM035-01	B01	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM035-01	B02	HmiRQP0168	MIMAT0000447	hsa-miR-134	
QM035-01	B03	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM035-01	B04	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM035-01	B05	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM035-01	B06	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM035-01	B07	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM035-01	B08	HmiRQP0510	MIMAT0001635	hsa-miR-452-5p	
QM035-01	B09	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM035-01	B10	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM035-01	B11	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM035-01	B12	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM035-01	C01	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM035-01	C02	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM035-01	C03	HmiRQP0476	MIMAT0000733	hsa-miR-379-5p	
QM035-01	C04	HmiRQP0447	MIMAT0004683	hsa-miR-362-3p	
QM035-01	C05	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM035-01	C06	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM035-01	C07	HmiRQP0479	MIMAT0000736	hsa-miR-381	
QM035-01	C08	HmiRQP0537	MIMAT0002817	hsa-miR-495	
QM035-01	C09	HmiRQP0369	MIMAT0002890	hsa-miR-299-5p	
QM035-01	C10	HmiRQP0446	MIMAT0000703	hsa-miR-361-5p	
QM035-01	C11	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM035-01	C12	HmiRQP0850	MIMAT0000094	hsa-miR-95	
QM035-01	D01	HmiRQP0448	MIMAT0000705	hsa-miR-362-5p	
QM035-01	D02	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM035-01	D03	HmiRQP0368	MIMAT0000687	hsa-miR-299-3p	
QM035-01	D04	HmiRQP0455	MIMAT0001621	hsa-miR-369-5p	
QM035-01	D05	HmiRQP0534	MIMAT0003161	hsa-miR-493-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM035-01	D06	HmiRQP0237	MIMAT0002821	hsa-miR-181d	
QM035-01	D07	HmiRQP0480	MIMAT0000737	hsa-miR-382-5p	
QM035-01	D08	HmiRQP0546	MIMAT0002872	hsa-miR-501-5p	
QM035-01	D09	HmiRQP0419	MIMAT0004700	hsa-miR-331-5p	
QM035-01	D10	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM035-01	D11	HmiRQP0445	MIMAT0004682	hsa-miR-361-3p	
QM035-01	D12	HmiRQP0484	MIMAT0001639	hsa-miR-409-3p	
QM035-01	E01	HmiRQP0522	MIMAT0004762	hsa-miR-486-3p	
QM035-01	E02	HmiRQP0420	MIMAT0000760	hsa-miR-331-3p	
QM035-01	E03	HmiRQP0501	MIMAT0002814	hsa-miR-432-5p	
QM035-01	E04	HmiRQP0198	MIMAT0004766	hsa-miR-146b-3p	
QM035-01	E05	HmiRQP0492	MIMAT0004748	hsa-miR-423-5p	
QM035-01	E06	HmiRQP0520	MIMAT0002175	hsa-miR-485-5p	
QM035-01	E07	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	
QM035-01	E08	HmiRQP0523	MIMAT0002177	hsa-miR-486-5p	
QM035-01	E09	HmiRQP0580	MIMAT0002863	hsa-miR-518a-3p	
QM035-01	E10	HmiRQP0287	MIMAT0000227	hsa-miR-197-3p	
QM035-01	E11	HmiRQP0835	MIMAT0003218	hsa-miR-92b-3p	
QM035-01	E12	HmiRQP0525	MIMAT0003180	hsa-miR-487b	
QM035-01	F01	HmiRQP0204	MIMAT0000243	hsa-miR-148a-3p	
QM035-01	F02	HmiRQP0467	MIMAT0000729	hsa-miR-376a-3p	
QM035-01	F03	HmiRQP0206	MIMAT0000759	hsa-miR-148b-3p	
QM035-01	F04	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM035-01	F05	HmiRQP0547	MIMAT0004774	hsa-miR-501-3p	
QM035-01	F06	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM035-01	F07	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM035-01	F08	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	hsa-miR-199b-3p
QM035-01	F09	HmiRQP0218	MIMAT0000453	hsa-miR-154-3p	
QM035-01	F10	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM035-01	F11	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM035-01	F12	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM035-01	G01	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM035-01	G02	HmiRQP0321	MIMAT0004564	hsa-miR-214-5p	
QM035-01	G03	HmiRQP0333	MIMAT0004495	hsa-miR-22-5p	
QM035-01	G04	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM035-01	G05	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM035-01	G06	HmiRQP0376	MIMAT0004673	hsa-miR-29c-5p	
QM035-01	G07	HmiRQP0422	MIMAT0004703	hsa-miR-335-3p	
QM035-01	G08	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM035-01	G09	HmiRQP0205	MIMAT0004699	hsa-miR-148b-5p	
QM035-01	G10	HmiRQP0279	MIMAT0004767	hsa-miR-193b-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM035-01	G11	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM035-01	G12	HmiRQP0340	MIMAT0004569	hsa-miR-222-5p	
QM035-01	H01	NC	NC	NC	
QM035-01	H02	NC	NC	NC	
QM035-01	H03	HmiRQP9001	NR_002752	HK1	
QM035-01	H04	HmiRQP9011	NR_002750	HK2	
QM035-01	H05	HmiRQP9021	NR_002745	HK3	
QM035-01	H06	HmiRQP9051	NR_002746	HK4	
QM035-01	H07	HmiRQP9061	NR_002744	HK5	
QM035-01	H08	HmiRQP9071	NR_002450	HK6	
QM035-01	H09	RT	RT	RT	
QM035-01	H10	RT	RT	RT	
QM035-01	H11	PCR	PCR	PCR	
QM035-01	H12	PCR	PCR	PCR	