

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM012-01	A01	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM012-01	A02	HmiRQP0029	MIMAT0000680	hsa-miR-106b-5p	
QM012-01	A03	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM012-01	A04	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM012-01	A05	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM012-01	A06	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM012-01	A07	HmiRQP0175	MIMAT0000429	hsa-miR-137	
QM012-01	A08	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM012-01	A09	HmiRQP0214	MIMAT0000439	hsa-miR-153	
QM012-01	A10	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM012-01	A11	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM012-01	A12	HmiRQP0258	MIMAT0000458	hsa-miR-190a	
QM012-01	B01	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM012-01	B02	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM012-01	B03	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM012-01	B04	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM012-01	B05	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM012-01	B06	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM012-01	B07	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM012-01	B08	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM012-01	B09	HmiRQP0317	MIMAT0000267	hsa-miR-210	
QM012-01	B10	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM012-01	B11	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM012-01	B12	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM012-01	C01	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM012-01	C02	HmiRQP0177	MIMAT0000430	hsa-miR-138-5p	
QM012-01	C03	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM012-01	C04	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM012-01	C05	HmiRQP0251	MIMAT0000262	hsa-miR-187-3p	
QM012-01	C06	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM012-01	C07	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM012-01	C08	HmiRQP0323	MIMAT0000273	hsa-miR-216a	
QM012-01	C09	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM012-01	C10	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM012-01	C11	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM012-01	C12	HmiRQP0364	MIMAT0000690	hsa-miR-296-5p	
QM012-01	D01	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM012-01	D02	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM012-01	D03	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM012-01	D04	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM012-01	D05	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM012-01	D06	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM012-01	D07	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM012-01	D08	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM012-01	D09	HmiRQP0325	MIMAT0000274	hsa-miR-217	
QM012-01	D10	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM012-01	D11	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM012-01	D12	HmiRQP0190	MIMAT0000436	hsa-miR-144-3p	
QM012-01	E01	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	
QM012-01	E02	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM012-01	E03	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM012-01	E04	HmiRQP0334	MIMAT0000277	hsa-miR-220a	
QM012-01	E05	HmiRQP0305	MIMAT0000264	hsa-miR-203	
QM012-01	E06	HmiRQP0362	MIMAT0000085	hsa-miR-28-5p	
QM012-01	E07	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM012-01	E08	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM012-01	E09	HmiRQP0125	MIMAT0000424	hsa-miR-128	
QM012-01	E10	HmiRQP0410	MIMAT0000755	hsa-miR-323a-3p	
QM012-01	E11	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM012-01	E12	HmiRQP0363	MIMAT0004502	hsa-miR-28-3p	
QM012-01	F01	HmiRQP0336	MIMAT0004915	hsa-miR-220c	
QM012-01	F02	HmiRQP0516	MIMAT0003150	hsa-miR-455-5p	
QM012-01	F03	HmiRQP0335	MIMAT0004908	hsa-miR-220b	
QM012-01	F04	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM012-01	F05	HmiRQP0515	MIMAT0004784	hsa-miR-455-3p	
QM012-01	F06	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM012-01	F07	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM012-01	F08	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM012-01	F09	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM012-01	F10	HmiRQP0412	MIMAT0000761	hsa-miR-324-5p	
QM012-01	F11	HmiRQP0419	MIMAT0004700	hsa-miR-331-5p	
QM012-01	F12	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM012-01	G01	HmiRQP0495	MIMAT0003393	hsa-miR-425-5p	
QM012-01	G02	HmiRQP0522	MIMAT0004762	hsa-miR-486-3p	
QM012-01	G03	HmiRQP0198	MIMAT0004766	hsa-miR-146b-3p	
QM012-01	G04	HmiRQP0414	MIMAT0000756	hsa-miR-326	
QM012-01	G05	HmiRQP0523	MIMAT0002177	hsa-miR-486-5p	
QM012-01	G06	HmiRQP0287	MIMAT0000227	hsa-miR-197-3p	
QM012-01	G07	HmiRQP0365	MIMAT0004679	hsa-miR-296-3p	
QM012-01	G08	HmiRQP0415	MIMAT0000752	hsa-miR-328	
QM012-01	G09	HmiRQP0409	MIMAT0004696	hsa-miR-323a-5p	
QM012-01	G10	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM012-01	G11	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM012-01	G12	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM012-01	H01	NC	NC	NC	
QM012-01	H02	NC	NC	NC	
QM012-01	H03	HmiRQP9001	NR_002752	HK1	
QM012-01	H04	HmiRQP9011	NR_002750	HK2	
QM012-01	H05	HmiRQP9021	NR_002745	HK3	
QM012-01	H06	HmiRQP9051	NR_002746	HK4	
QM012-01	H07	HmiRQP9061	NR_002744	HK5	
QM012-01	H08	HmiRQP9071	NR_002450	HK6	
QM012-01	H09	RT	RT	RT	
QM012-01	H10	RT	RT	RT	
QM012-01	H11	PCR	PCR	PCR	
QM012-01	H12	PCR	PCR	PCR	