

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM046_01	A01	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM046_01	A02	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM046_01	A03	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM046_01	A04	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM046_01	A05	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM046_01	A06	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM046_01	A07	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM046_01	A08	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM046_01	A09	HmiRQP0029	MIMAT0000680	hsa-miR-106b-5p	
QM046_01	A10	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM046_01	A11	HmiRQP0031	MIMAT0004555	hsa-miR-10a	
QM046_01	A12	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM046_01	B01	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM046_01	B02	HmiRQP3286	MIMAT0022713	hsa-miR-1185-2-3p	
QM046_01	B03	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM046_01	B04	HmiRQP0078	MIMAT0005898	hsa-miR-1246	
QM046_01	B05	HmiRQP0087	MIMAT0005906	hsa-miR-1255a	
QM046_01	B06	HmiRQP0093	MIMAT0004602	hsa-miR-125a	
QM046_01	B07	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM046_01	B08	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM046_01	B09	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM046_01	B10	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	
QM046_01	B11	HmiRQP0134	MIMAT0005880	hsa-miR-1290	
QM046_01	B12	HmiRQP0152	MIMAT0005893	hsa-miR-1305	
QM046_01	C01	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM046_01	C02	HmiRQP0160	MIMAT0004594	hsa-miR-132	
QM046_01	C03	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM046_01	C04	HmiRQP0173	MIMAT0000448	hsa-miR-136-5p	
QM046_01	C05	HmiRQP0175	MIMAT0000429	hsa-miR-137	
QM046_01	C06	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM046_01	C07	HmiRQP0186	MIMAT0000434	hsa-miR-142-3p	
QM046_01	C08	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM046_01	C09	HmiRQP0190	MIMAT0000436	hsa-miR-144-3p	
QM046_01	C10	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM046_01	C11	HmiRQP0195	MIMAT0004608	hsa-miR-146a	
QM046_01	C12	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM046_01	D01	HmiRQP0204	MIMAT0000243	hsa-miR-148a-3p	
QM046_01	D02	HmiRQP0206	MIMAT0000759	hsa-miR-148b-3p	
QM046_01	D03	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM046_01	D04	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM046_01	D05	HmiRQP0211	MIMAT0000757	hsa-miR-151a-3p	
QM046_01	D06	HmiRQP0219	MIMAT0000452	hsa-miR-154-5p	
QM046_01	D07	HmiRQP0222	MIMAT0004488	hsa-miR-15a	
QM046_01	D08	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM046_01	D09	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM046_01	D10	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM046_01	D11	HmiRQP0233	MIMAT0004558	hsa-miR-181a-2-3p	
QM046_01	D12	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM046_01	E01	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM046_01	E02	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM046_01	E03	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM046_01	E04	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM046_01	E05	HmiRQP0251	MIMAT0000262	hsa-miR-187-3p	
QM046_01	E06	HmiRQP0255	MIMAT0000072	hsa-miR-18a-5p	
QM046_01	E07	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM046_01	E08	HmiRQP0277	MIMAT0004614	hsa-miR-193a-5p	
QM046_01	E09	HmiRQP0279	MIMAT0004767	hsa-miR-193b-5p	
QM046_01	E10	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM046_01	E11	HmiRQP0286	MIMAT0001080	hsa-miR-196b-5p	
QM046_01	E12	HmiRQP3038	MIMAT0022691	hsa-miR-197-5p	
QM046_01	F01	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM046_01	F02	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM046_01	F03	HmiRQP0289	MIMAT0004563	hsa-miR-199b-3p	
QM046_01	F04	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM046_01	F05	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM046_01	F06	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM046_01	F07	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	
QM046_01	F08	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM046_01	F09	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM046_01	F10	HmiRQP0314	MIMAT0001413	hsa-miR-20b-5p	
QM046_01	F11	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM046_01	F12	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM046_01	G01	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM046_01	G02	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM046_01	G03	HmiRQP0339	MIMAT0000279	hsa-miR-222	
QM046_01	G04	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM046_01	G05	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM046_01	G06	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM046_01	G07	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM046_01	G08	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM046_01	G09	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM046_01	G10	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM046_01	G11	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM046_01	G12	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM046_01	H01	NC	NC	NC	
QM046_01	H02	NC	NC	NC	
QM046_01	H03	HKG1	HKG1	HKG1	
QM046_01	H04	HKG2	HKG2	HKG2	
QM046_01	H05	HKG3	HKG3	HKG3	
QM046_01	H06	HKG4	HKG4	HKG4	
QM046_01	H07	HKG5	HKG5	HKG5	
QM046_01	H08	HKG6	HKG6	HKG6	
QM046_01	H09	RT	RT	RT	
QM046_01	H10	RT	RT	RT	
QM046_01	H11	PCR	PCR	PCR	
QM046_01	H12	PCR	PCR	PCR	
QM046_02	A01	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM046_02	A02	HmiRQP0362	MIMAT0000085	hsa-miR-28-5p	
QM046_02	A03	HmiRQP0364	MIMAT0000690	hsa-miR-296-5p	
QM046_02	A04	HmiRQP0369	MIMAT0002890	hsa-miR-299-5p	
QM046_02	A05	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM046_02	A06	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM046_02	A07	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM046_02	A08	HmiRQP0378	MIMAT0000688	hsa-miR-301a-3p	
QM046_02	A09	HmiRQP1518	MIMAT0015066	hsa-miR-3065-5p	
QM046_02	A10	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM046_02	A11	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM046_02	A12	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM046_02	B01	HmiRQP1523	MIMAT0015072	hsa-miR-320e	
QM046_02	B02	HmiRQP0410	MIMAT0000755	hsa-miR-323a-3p	
QM046_02	B03	HmiRQP0404	MIMAT0000090	hsa-miR-32-5p	
QM046_02	B04	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM046_02	B05	HmiRQP0423	MIMAT0000754	hsa-miR-337-3p	
QM046_02	B06	HmiRQP0424	MIMAT0004695	hsa-miR-337-5p	
QM046_02	B07	HmiRQP0436	MIMAT0000753	hsa-miR-342-3p	
QM046_02	B08	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM046_02	B09	HmiRQP1907	MIMAT0017991	hsa-miR-3613-3p	
QM046_02	B10	HmiRQP0446	MIMAT0000703	hsa-miR-361-5p	
QM046_02	B11	HmiRQP0448	MIMAT0000705	hsa-miR-362-5p	
QM046_02	B12	HmiRQP0449	MIMAT0000707	hsa-miR-363-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM046_02	C01	HmiRQP0451	MIMAT0000710	hsa-miR-365a-3p	
QM046_02	C02	HmiRQP1944	MIMAT0018087	hsa-miR-3665	
QM046_02	C03	HmiRQP0454	MIMAT0000721	hsa-miR-369-3p	
QM046_02	C04	HmiRQP0463	MIMAT0000727	hsa-miR-374a-5p	
QM046_02	C05	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM046_02	C06	HmiRQP0467	MIMAT0000729	hsa-miR-376a-3p	
QM046_02	C07	HmiRQP0472	MIMAT0000730	hsa-miR-377-3p	
QM046_02	C08	HmiRQP2046	MIMAT0018937	hsa-miR-378g	
QM046_02	C09	HmiRQP2182	MIMAT0019074	hsa-miR-378i	
QM046_02	C10	HmiRQP0476	MIMAT0000733	hsa-miR-379-5p	
QM046_02	C11	HmiRQP0480	MIMAT0000737	hsa-miR-382-5p	
QM046_02	C12	HmiRQP0484	MIMAT0001639	hsa-miR-409-3p	
QM046_02	D01	HmiRQP0487	MIMAT0004813	hsa-miR-411-3p	
QM046_02	D02	HmiRQP0486	MIMAT0003329	hsa-miR-411-5p	
QM046_02	D03	HmiRQP0492	MIMAT0004748	hsa-miR-423-5p	
QM046_02	D04	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM046_02	D05	HmiRQP1612	MIMAT0016916	hsa-miR-4286	
QM046_02	D06	HmiRQP0501	MIMAT0002814	hsa-miR-432-5p	
QM046_02	D07	HmiRQP2084	MIMAT0018976	hsa-miR-4454	
QM046_02	D08	HmiRQP2100	MIMAT0018993	hsa-miR-4466	
QM046_02	D09	HmiRQP0503	MIMAT0001532	hsa-miR-448	
QM046_02	D10	HmiRQP2128	MIMAT0019022	hsa-miR-4488	
QM046_02	D11	HmiRQP0506	MIMAT0001545	hsa-miR-450a-5p	
QM046_02	D12	HmiRQP2159	MIMAT0019053	hsa-miR-4516	
QM046_02	E01	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM046_02	E02	HmiRQP0516	MIMAT0003150	hsa-miR-455-5p	
QM046_02	E03	HmiRQP2241	MIMAT0019709	hsa-miR-4647	
QM046_02	E04	HmiRQP2495	MIMAT0019964	hsa-miR-4792	
QM046_02	E05	HmiRQP0540	MIMAT0002820	hsa-miR-497-5p	
QM046_02	E06	HmiRQP0543	MIMAT0002870	hsa-miR-499a-5p	
QM046_02	E07	HmiRQP1537	MIMAT0015087	hsa-miR-514b-5p	
QM046_02	E08	HmiRQP0610	MIMAT0002838	hsa-miR-525-5p	
QM046_02	E09	HmiRQP0616	MIMAT0004780	hsa-miR-532-3p	
QM046_02	E10	HmiRQP0622	MIMAT0004954	hsa-miR-543	
QM046_02	E11	HmiRQP0627	MIMAT0004803	hsa-miR-548a-5p	
QM046_02	E12	HmiRQP2631	MIMAT0022266	hsa-miR-548ar-3p	
QM046_02	F01	HmiRQP2015	MIMAT0018354	hsa-miR-548y	
QM046_02	F02	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM046_02	F03	HmiRQP0674	MIMAT0004795	hsa-miR-574-5p	
QM046_02	F04	HmiRQP0675	MIMAT0003240	hsa-miR-575	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM046_02	F05	HmiRQP0684	MIMAT0003247	hsa-miR-582-5p	
QM046_02	F06	HmiRQP2775	MIMAT0023714	hsa-miR-6089	
QM046_02	F07	HmiRQP0716	MIMAT0003280	hsa-miR-612	
QM046_02	F08	HmiRQP0720	MIMAT0003283	hsa-miR-615-3p	
QM046_02	F09	HmiRQP0734	MIMAT0003295	hsa-miR-626	
QM046_02	F10	HmiRQP0740	MIMAT0003299	hsa-miR-630	
QM046_02	F11	HmiRQP0760	MIMAT0003320	hsa-miR-650	
QM046_02	F12	HmiRQP0769	MIMAT0003336	hsa-miR-658	
QM046_02	G01	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM046_02	G02	HmiRQP0796	MIMAT0003882	hsa-miR-767-5p	
QM046_02	G03	HmiRQP4563	MIMAT0031178	hsa-miR-7975	
QM046_02	G04	HmiRQP0802	MIMAT0004185	hsa-miR-802	
QM046_02	G05	HmiRQP0817	MIMAT0004916	hsa-miR-888-5p	
QM046_02	G06	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM046_02	G07	HmiRQP0853	MIMAT0000096	hsa-miR-98-5p	
QM046_02	G08	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM046_02	G09	HmiRQP0856	MIMAT0004678	hsa-miR-99b	
QM046_02	G10	HmiRQP0857	MIMAT0000689	hsa-miR-99b-5p	
QM046_02	H01	NC	NC	NC	
QM046_02	H02	NC	NC	NC	
QM046_02	H03	HKG1	HKG1	HKG1	
QM046_02	H04	HKG2	HKG2	HKG2	
QM046_02	H05	HKG3	HKG3	HKG3	
QM046_02	H06	HKG4	HKG4	HKG4	
QM046_02	H07	HKG5	HKG5	HKG5	
QM046_02	H08	HKG6	HKG6	HKG6	
QM046_02	H09	RT	RT	RT	
QM046_02	H10	RT	RT	RT	
QM046_02	H11	PCR	PCR	PCR	
QM046_02	H12	PCR	PCR	PCR	