

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM017-01	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1	
QM017-01	A02	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM017-01	A03	HmiRQP0029	MIMAT0000680	hsa-miR-106b-5p	
QM017-01	A04	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM017-01	A05	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM017-01	A06	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM017-01	A07	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM017-01	A08	HmiRQP0214	MIMAT0000439	hsa-miR-153	
QM017-01	A09	HmiRQP0219	MIMAT0000452	hsa-miR-154-5p	
QM017-01	A10	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM017-01	A11	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM017-01	A12	HmiRQP0258	MIMAT0000458	hsa-miR-190a	
QM017-01	B01	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM017-01	B02	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM017-01	B03	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM017-01	B04	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM017-01	B05	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM017-01	B06	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM017-01	B07	HmiRQP0317	MIMAT0000267	hsa-miR-210	
QM017-01	B08	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM017-01	B09	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM017-01	B10	HmiRQP0459	MIMAT0000724	hsa-miR-372	
QM017-01	B11	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM017-01	B12	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM017-01	C01	HmiRQP0168	MIMAT0000447	hsa-miR-134	
QM017-01	C02	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM017-01	C03	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM017-01	C04	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM017-01	C05	HmiRQP0252	MIMAT0000457	hsa-miR-188-5p	
QM017-01	C06	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM017-01	C07	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM017-01	C08	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM017-01	C09	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM017-01	C10	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM017-01	C11	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM017-01	C12	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM017-01	D01	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM017-01	D02	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM017-01	D03	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM017-01	D04	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM017-01	D05	HmiRQP0325	MIMAT0000274	hsa-miR-217	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM017-01	D06	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM017-01	D07	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM017-01	D08	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	
QM017-01	D09	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM017-01	D10	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM017-01	D11	HmiRQP0434	MIMAT0004692	hsa-miR-340-5p	
QM017-01	D12	HmiRQP0413	MIMAT0000771	hsa-miR-325	
QM017-01	E01	HmiRQP0334	MIMAT0000277	hsa-miR-220a	
QM017-01	E02	HmiRQP0305	MIMAT0000264	hsa-miR-203	
QM017-01	E03	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM017-01	E04	HmiRQP0212	MIMAT0004697	hsa-miR-151a-5p	
QM017-01	E05	HmiRQP0125	MIMAT0000424	hsa-miR-128	
QM017-01	E06	HmiRQP0211	MIMAT0000757	hsa-miR-151a-3p	
QM017-01	E07	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM017-01	E08	HmiRQP0435	MIMAT0004694	hsa-miR-342-5p	
QM017-01	E09	HmiRQP0336	MIMAT0004915	hsa-miR-220c	
QM017-01	E10	HmiRQP0199	MIMAT0000251	hsa-miR-147a	
QM017-01	E11	HmiRQP0452	MIMAT0000719	hsa-miR-367-3p	
QM017-01	E12	HmiRQP0430	MIMAT0000091	hsa-miR-33a-5p	
QM017-01	F01	HmiRQP0335	MIMAT0004908	hsa-miR-220b	
QM017-01	F02	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM017-01	F03	HmiRQP0783	MIMAT0004926	hsa-miR-708-5p	
QM017-01	F04	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM017-01	F05	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM017-01	F06	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM017-01	F07	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM017-01	F08	HmiRQP0412	MIMAT0000761	hsa-miR-324-5p	
QM017-01	F09	HmiRQP0419	MIMAT0004700	hsa-miR-331-5p	
QM017-01	F10	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM017-01	F11	HmiRQP0420	MIMAT0000760	hsa-miR-331-3p	
QM017-01	F12	HmiRQP0436	MIMAT0000753	hsa-miR-342-3p	
QM017-01	G01	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM017-01	G02	HmiRQP0414	MIMAT0000756	hsa-miR-326	
QM017-01	G03	HmiRQP0253	MIMAT0004613	hsa-miR-188-3p	
QM017-01	G04	HmiRQP0415	MIMAT0000752	hsa-miR-328	
QM017-01	G05	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM017-01	G06	HmiRQP0463	MIMAT0000727	hsa-miR-374a-5p	
QM017-01	G07	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM017-01	G08	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM017-01	G09	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM017-01	G10	HmiRQP0853	MIMAT0000096	hsa-miR-98	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM017-01	G11	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM017-01	G12	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM017-01	H01	NC	NC	NC	
QM017-01	H02	NC	NC	NC	
QM017-01	H03	HmiRQP9001	NR_002752	HK1	
QM017-01	H04	HmiRQP9011	NR_002750	HK2	
QM017-01	H05	HmiRQP9021	NR_002745	HK3	
QM017-01	H06	HmiRQP9051	NR_002746	HK4	
QM017-01	H07	HmiRQP9061	NR_002744	HK5	
QM017-01	H08	HmiRQP9071	NR_002450	HK6	
QM017-01	H09	RT	RT	RT	
QM017-01	H10	RT	RT	RT	
QM017-01	H11	PCR	PCR	PCR	
QM017-01	H12	PCR	PCR	PCR	
QM017-02	A01	HmiRQP0408	MIMAT0006764	hsa-miR-320d	
QM017-02	A02	HmiRQP0387	MIMAT0000718	hsa-miR-302d-3p	
QM017-02	A03	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM017-02	A04	HmiRQP0314	MIMAT0001413	hsa-miR-20b-5p	
QM017-02	A05	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM017-02	A06	HmiRQP0407	MIMAT0005793	hsa-miR-320c	
QM017-02	A07	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM017-02	A08	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM017-02	A09	HmiRQP0159	MIMAT0000691	hsa-miR-130b-3p	
QM017-02	A10	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM017-02	A11	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM017-02	A12	HmiRQP0006	MIMAT0000064	hsa-let-7c	
QM017-02	B01	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM017-02	B02	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM017-02	B03	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM017-02	B04	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM017-02	B05	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM017-02	B06	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM017-02	B07	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM017-02	B08	HmiRQP0286	MIMAT0001080	hsa-miR-196b-5p	
QM017-02	B09	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM017-02	B10	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM017-02	B11	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM017-02	B12	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM017-02	C01	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM017-02	C02	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM017-02	C03	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM017-02	C04	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM017-02	C05	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM017-02	C06	HmiRQP0405	MIMAT0000510	hsa-miR-320a	
QM017-02	C07	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM017-02	C08	HmiRQP0462	MIMAT0004688	hsa-miR-374a-3p	
QM017-02	C09	HmiRQP0236	MIMAT0004559	hsa-miR-181c-3p	
QM017-02	C10	HmiRQP0275	MIMAT0004543	hsa-miR-192-3p	
QM017-02	C11	HmiRQP0233	MIMAT0004558	hsa-miR-181a-2-3p	
QM017-02	C12	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM017-02	D01	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM017-02	D02	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM017-02	D03	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM017-02	D04	HmiRQP0005	MIMAT0004483	hsa-let-7c*	
QM017-02	D05	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM017-02	D06	HmiRQP0009	MIMAT0004485	hsa-let-7e-3p	
QM017-02	D07	HmiRQP0011	MIMAT0004486	hsa-let-7f-1-3p	
QM017-02	D08	HmiRQP0013	MIMAT0004487	hsa-let-7f-2-3p	
QM017-02	D09	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM017-02	D10	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM017-02	D11	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM017-02	D12	HmiRQP0033	MIMAT0004556	hsa-miR-10b-3p	
QM017-02	E01	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM017-02	E02	HmiRQP0158	MIMAT0004680	hsa-miR-130b-5p	
QM017-02	E03	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM017-02	E04	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM017-02	E05	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM017-02	E06	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM017-02	E07	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM017-02	E08	HmiRQP0333	MIMAT0004495	hsa-miR-22-5p	
QM017-02	E09	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM017-02	E10	HmiRQP0345	MIMAT0004496	hsa-miR-23a-5p	
QM017-02	E11	HmiRQP0347	MIMAT0004587	hsa-miR-23b-5p	
QM017-02	E12	HmiRQP0348	MIMAT0000079	hsa-miR-24-1-5p	
QM017-02	F01	HmiRQP0360	MIMAT0004588	hsa-miR-27b-5p	
QM017-02	F02	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM017-02	F03	HmiRQP0372	MIMAT0004514	hsa-miR-29b-1-5p	
QM017-02	F04	HmiRQP0374	MIMAT0004515	hsa-miR-29b-2-5p	
QM017-02	F05	HmiRQP0376	MIMAT0004673	hsa-miR-29c-5p	
QM017-02	F06	HmiRQP0397	MIMAT0004551	hsa-miR-30d-3p	
QM017-02	F07	HmiRQP0422	MIMAT0004703	hsa-miR-335-3p	
QM017-02	F08	HmiRQP0429	MIMAT0004506	hsa-miR-33a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM017-02	F09	HmiRQP0442	MIMAT0000685	hsa-miR-34b-5p	
QM017-02	F10	HmiRQP0453	MIMAT0004686	hsa-miR-367-5p	
QM017-02	F11	HmiRQP0465	MIMAT0004956	hsa-miR-374b-3p	
QM017-02	F12	HmiRQP0493	MIMAT0004749	hsa-miR-424-3p	
QM017-02	G01	HmiRQP0784	MIMAT0004927	hsa-miR-708-3p	
QM017-02	G02	HmiRQP1309	MIMAT0010195	hsa-let-7a-2-3p	
QM017-02	G03	HmiRQP1523	MIMAT0015072	hsa-miR-320e	
QM017-02	G04	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM017-02	G05	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM017-02	G06	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM017-02	G07	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM017-02	G08	HmiRQP0224	MIMAT0004586	hsa-miR-15b-3p	
QM017-02	G09	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM017-02	G10	HmiRQP0228	MIMAT0004518	hsa-miR-16-2-3p	
QM017-02	G11	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM017-02	G12	HmiRQP0243	MIMAT0004560	hsa-miR-183-3p	
QM017-02	H01	NC	NC	NC	
QM017-02	H02	NC	NC	NC	
QM017-02	H03	HmiRQP9001	NR_002752	HK1	
QM017-02	H04	HmiRQP9011	NR_002750	HK2	
QM017-02	H05	HmiRQP9021	NR_002745	HK3	
QM017-02	H06	HmiRQP9051	NR_002746	HK4	
QM017-02	H07	HmiRQP9061	NR_002744	HK5	
QM017-02	H08	HmiRQP9071	NR_002450	HK6	
QM017-02	H09	RT	RT	RT	
QM017-02	H10	RT	RT	RT	
QM017-02	H11	PCR	PCR	PCR	
QM017-02	H12	PCR	PCR	PCR	