

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM009-01	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1	
QM009-01	A02	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM009-01	A03	HmiRQP0029	MIMAT0000680	hsa-miR-106b-5p	
QM009-01	A04	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM009-01	A05	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM009-01	A06	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM009-01	A07	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM009-01	A08	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM009-01	A09	HmiRQP0175	MIMAT0000429	hsa-miR-137	
QM009-01	A10	HmiRQP0180	MIMAT0000250	hsa-miR-139-5p	
QM009-01	A11	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM009-01	A12	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM009-01	B01	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM009-01	B02	HmiRQP0330	MIMAT0000276	hsa-miR-219-5p	
QM009-01	B03	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM009-01	B04	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM009-01	B05	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM009-01	B06	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM009-01	B07	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM009-01	B08	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM009-01	B09	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM009-01	B10	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM009-01	B11	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM009-01	B12	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM009-01	C01	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM009-01	C02	HmiRQP0186	MIMAT0000434	hsa-miR-142-3p	
QM009-01	C03	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM009-01	C04	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM009-01	C05	HmiRQP0213	MIMAT0000438	hsa-miR-152	
QM009-01	C06	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM009-01	C07	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM009-01	C08	HmiRQP0280	MIMAT0000460	hsa-miR-194-5p	
QM009-01	C09	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM009-01	C10	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM009-01	C11	HmiRQP0323	MIMAT0000273	hsa-miR-216a	
QM009-01	C12	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM009-01	D01	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM009-01	D02	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM009-01	D03	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM009-01	D04	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM009-01	D05	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM009-01	D06	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM009-01	D07	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM009-01	D08	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM009-01	D09	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM009-01	D10	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM009-01	D11	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM009-01	D12	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM009-01	E01	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM009-01	E02	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM009-01	E03	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM009-01	E04	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM009-01	E05	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM009-01	E06	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM009-01	E07	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM009-01	E08	HmiRQP0305	MIMAT0000264	hsa-miR-203	
QM009-01	E09	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM009-01	E10	HmiRQP0212	MIMAT0004697	hsa-miR-151a-5p	
QM009-01	E11	HmiRQP0125	MIMAT0000424	hsa-miR-128	
QM009-01	E12	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM009-01	F01	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM009-01	F02	HmiRQP0425	MIMAT0004701	hsa-miR-338-5p	
QM009-01	F03	HmiRQP0324	MIMAT0004959	hsa-miR-216b	
QM009-01	F04	HmiRQP0545	MIMAT0004773	hsa-miR-500a-5p	
QM009-01	F05	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM009-01	F06	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM009-01	F07	HmiRQP0139	MIMAT0004605	hsa-miR-129-2-3p	
QM009-01	F08	HmiRQP0412	MIMAT0000761	hsa-miR-324-5p	
QM009-01	F09	HmiRQP0426	MIMAT0000763	hsa-miR-338-3p	
QM009-01	F10	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM009-01	F11	HmiRQP0329	MIMAT0004567	hsa-miR-219-1-3p	
QM009-01	F12	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM009-01	G01	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	
QM009-01	G02	HmiRQP0415	MIMAT0000752	hsa-miR-328	
QM009-01	G03	HmiRQP0179	MIMAT0004552	hsa-miR-139-3p	
QM009-01	G04	HmiRQP0456	MIMAT0000722	hsa-miR-370	
QM009-01	G05	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM009-01	G06	HmiRQP0463	MIMAT0000727	hsa-miR-374a-5p	
QM009-01	G07	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM009-01	G08	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM009-01	G09	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM009-01	G10	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM009-01	G11	HmiRQP0204	MIMAT0000243	hsa-miR-148a-3p	
QM009-01	G12	HmiRQP0853	MIMAT0000096	hsa-miR-98	
QM009-01	H01	NC	NC	NC	
QM009-01	H02	NC	NC	NC	
QM009-01	H03	HmiRQP9001	NR_002752	HK1	
QM009-01	H04	HmiRQP9011	NR_002750	HK2	
QM009-01	H05	HmiRQP9021	NR_002745	HK3	
QM009-01	H06	HmiRQP9051	NR_002746	HK4	
QM009-01	H07	HmiRQP9061	NR_002744	HK5	
QM009-01	H08	HmiRQP9071	NR_002450	HK6	
QM009-01	H09	RT	RT	RT	
QM009-01	H10	RT	RT	RT	
QM009-01	H11	PCR	PCR	PCR	
QM009-01	H12	PCR	PCR	PCR	
QM009-02	A01	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM009-02	A02	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM009-02	A03	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM009-02	A04	HmiRQP0206	MIMAT0000759	hsa-miR-148b-3p	
QM009-02	A05	HmiRQP0322	MIMAT0000272	hsa-miR-215	
QM009-02	A06	HmiRQP0408	MIMAT0006764	hsa-miR-320d	
QM009-02	A07	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM009-02	A08	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM009-02	A09	HmiRQP0378	MIMAT0000688	hsa-miR-301a-3p	
QM009-02	A10	HmiRQP0407	MIMAT0005793	hsa-miR-320c	
QM009-02	A11	HmiRQP0379	MIMAT0004958	hsa-miR-301b	
QM009-02	A12	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM009-02	B01	HmiRQP0159	MIMAT0000691	hsa-miR-130b-3p	
QM009-02	B02	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM009-02	B03	HmiRQP0006	MIMAT0000064	hsa-let-7c	
QM009-02	B04	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM009-02	B05	HmiRQP0255	MIMAT0000072	hsa-miR-18a-5p	
QM009-02	B06	HmiRQP0026	MIMAT0000103	hsa-miR-106a-5p	
QM009-02	B07	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM009-02	B08	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM009-02	B09	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM009-02	B10	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM009-02	B11	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM009-02	B12	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM009-02	C01	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM009-02	C02	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM009-02	C03	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM009-02	C04	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM009-02	C05	HmiRQP0405	MIMAT0000510	hsa-miR-320a	
QM009-02	C06	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM009-02	C07	HmiRQP0166	MIMAT0000427	hsa-miR-133a	
QM009-02	C08	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM009-02	C09	HmiRQP0462	MIMAT0004688	hsa-miR-374a-3p	
QM009-02	C10	HmiRQP0236	MIMAT0004559	hsa-miR-181c-3p	
QM009-02	C11	HmiRQP0275	MIMAT0004543	hsa-miR-192-3p	
QM009-02	C12	HmiRQP0233	MIMAT0004558	hsa-miR-181a-2-3p	
QM009-02	D01	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM009-02	D02	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM009-02	D03	HmiRQP0299	MIMAT0004571	hsa-miR-200b-5p	
QM009-02	D04	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM009-02	D05	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM009-02	D06	HmiRQP0005	MIMAT0004483	hsa-let-7c*	
QM009-02	D07	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM009-02	D08	HmiRQP0009	MIMAT0004485	hsa-let-7e-3p	
QM009-02	D09	HmiRQP0011	MIMAT0004486	hsa-let-7f-1-3p	
QM009-02	D10	HmiRQP0013	MIMAT0004487	hsa-let-7f-2-3p	
QM009-02	D11	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM009-02	D12	HmiRQP0019	MIMAT0004512	hsa-miR-100-3p	
QM009-02	E01	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM009-02	E02	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM009-02	E03	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM009-02	E04	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM009-02	E05	HmiRQP0136	MIMAT0004548	hsa-miR-129-1-3p	
QM009-02	E06	HmiRQP0158	MIMAT0004680	hsa-miR-130b-5p	
QM009-02	E07	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM009-02	E08	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM009-02	E09	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM009-02	E10	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM009-02	E11	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM009-02	E12	HmiRQP0249	MIMAT0004612	hsa-miR-186-3p	
QM009-02	F01	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM009-02	F02	HmiRQP0321	MIMAT0004564	hsa-miR-214-5p	
QM009-02	F03	HmiRQP0333	MIMAT0004495	hsa-miR-22-5p	
QM009-02	F04	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM009-02	F05	HmiRQP0345	MIMAT0004496	hsa-miR-23a-5p	
QM009-02	F06	HmiRQP0347	MIMAT0004587	hsa-miR-23b-5p	
QM009-02	F07	HmiRQP0350	MIMAT0004497	hsa-miR-24-2-5p	
QM009-02	F08	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM009-02	F09	HmiRQP0353	MIMAT0004499	hsa-miR-26a-1-3p	
QM009-02	F10	HmiRQP0355	MIMAT0004681	hsa-miR-26a-2-3p	
QM009-02	F11	HmiRQP0394	MIMAT0004674	hsa-miR-30c-1-3p	
QM009-02	F12	HmiRQP0402	MIMAT0004504	hsa-miR-31-3p	
QM009-02	G01	HmiRQP0465	MIMAT0004956	hsa-miR-374b-3p	
QM009-02	G02	HmiRQP0544	MIMAT0002871	hsa-miR-500a-3p	
QM009-02	G03	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM009-02	G04	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM009-02	G05	HmiRQP0855	MIMAT0004511	hsa-miR-99a-3p	
QM009-02	G06	HmiRQP1295	MIMAT0009198	hsa-miR-224-3p	
QM009-02	G07	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM009-02	G08	HmiRQP1309	MIMAT0010195	hsa-let-7a-2-3p	
QM009-02	G09	HmiRQP1523	MIMAT0015072	hsa-miR-320e	
QM009-02	G10	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM009-02	G11	HmiRQP0097	MIMAT0004603	hsa-miR-125b-2-3p	
QM009-02	G12	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM009-02	H01	NC	NC	NC	
QM009-02	H02	NC	NC	NC	
QM009-02	H03	HmiRQP9001	NR_002752	HK1	
QM009-02	H04	HmiRQP9011	NR_002750	HK2	
QM009-02	H05	HmiRQP9021	NR_002745	HK3	
QM009-02	H06	HmiRQP9051	NR_002746	HK4	
QM009-02	H07	HmiRQP9061	NR_002744	HK5	
QM009-02	H08	HmiRQP9071	NR_002450	HK6	
QM009-02	H09	RT	RT	RT	
QM009-02	H10	RT	RT	RT	
QM009-02	H11	PCR	PCR	PCR	
QM009-02	H12	PCR	PCR	PCR	