

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM005-01	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1	
QM005-01	A02	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM005-01	A03	HmiRQP0025	MIMAT0000102	hsa-miR-105-5p	
QM005-01	A04	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM005-01	A05	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM005-01	A06	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM005-01	A07	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM005-01	A08	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM005-01	A09	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM005-01	A10	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM005-01	A11	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM005-01	A12	HmiRQP0258	MIMAT0000458	hsa-miR-190a	
QM005-01	B01	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM005-01	B02	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM005-01	B03	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM005-01	B04	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM005-01	B05	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM005-01	B06	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM005-01	B07	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM005-01	B08	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM005-01	B09	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM005-01	B10	HmiRQP0317	MIMAT0000267	hsa-miR-210	
QM005-01	B11	HmiRQP0319	MIMAT0000269	hsa-miR-212-3p	
QM005-01	B12	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM005-01	C01	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM005-01	C02	HmiRQP0173	MIMAT0000448	hsa-miR-136-5p	
QM005-01	C03	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM005-01	C04	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM005-01	C05	HmiRQP0213	MIMAT0000438	hsa-miR-152	
QM005-01	C06	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM005-01	C07	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM005-01	C08	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM005-01	C09	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM005-01	C10	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM005-01	C11	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM005-01	C12	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM005-01	D01	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM005-01	D02	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM005-01	D03	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM005-01	D04	HmiRQP0461	MIMAT0000726	hsa-miR-373-3p	
QM005-01	D05	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM005-01	D06	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM005-01	D07	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM005-01	D08	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM005-01	D09	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM005-01	D10	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM005-01	D11	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM005-01	D12	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM005-01	E01	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM005-01	E02	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM005-01	E03	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM005-01	E04	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM005-01	E05	HmiRQP0540	MIMAT0002820	hsa-miR-497-5p	
QM005-01	E06	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM005-01	E07	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM005-01	E08	HmiRQP0434	MIMAT0004692	hsa-miR-340-5p	
QM005-01	E09	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM005-01	E10	HmiRQP0305	MIMAT0000264	hsa-miR-203	
QM005-01	E11	HmiRQP0451	MIMAT0000710	hsa-miR-365a-3p	
QM005-01	E12	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM005-01	F01	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM005-01	F02	HmiRQP0574	MIMAT0006778	hsa-miR-516a-3p	
QM005-01	F03	HmiRQP0537	MIMAT0002817	hsa-miR-495	
QM005-01	F04	HmiRQP0125	MIMAT0000424	hsa-miR-128	
QM005-01	F05	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM005-01	F06	HmiRQP0561	MIMAT0002882	hsa-miR-510	
QM005-01	F07	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM005-01	F08	HmiRQP0717	MIMAT0003281	hsa-miR-613	
QM005-01	F09	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM005-01	F10	HmiRQP0528	MIMAT0002805	hsa-miR-489	
QM005-01	F11	HmiRQP0303	MIMAT0002811	hsa-miR-202-3p	
QM005-01	F12	HmiRQP0237	MIMAT0002821	hsa-miR-181d	
QM005-01	G01	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM005-01	G02	HmiRQP0198	MIMAT0004766	hsa-miR-146b-3p	
QM005-01	G03	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM005-01	G04	HmiRQP0520	MIMAT0002175	hsa-miR-485-5p	
QM005-01	G05	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	
QM005-01	G06	HmiRQP0415	MIMAT0000752	hsa-miR-328	
QM005-01	G07	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM005-01	G08	HmiRQP0774	MIMAT0003326	hsa-miR-663a	
QM005-01	G09	HmiRQP0772	MIMAT0003324	hsa-miR-661	
QM005-01	G10	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM005-01	G11	HmiRQP0017	MIMAT0000415	hsa-let-7i-5p	
QM005-01	G12	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM005-01	H01	NC	NC	NC	
QM005-01	H02	NC	NC	NC	
QM005-01	H03	HmiRQP9001	NR_002752	HK1	
QM005-01	H04	HmiRQP9011	NR_002750	HK2	
QM005-01	H05	HmiRQP9021	NR_002745	HK3	
QM005-01	H06	HmiRQP9051	NR_002746	HK4	
QM005-01	H07	HmiRQP9061	NR_002744	HK5	
QM005-01	H08	HmiRQP9071	NR_002450	HK6	
QM005-01	H09	RT	RT	RT	
QM005-01	H10	RT	RT	RT	
QM005-01	H11	PCR	PCR	PCR	
QM005-01	H12	PCR	PCR	PCR	
QM005-02	A01	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	
QM005-02	A02	HmiRQP0204	MIMAT0000243	hsa-miR-148a-3p	
QM005-02	A03	HmiRQP0853	MIMAT0000096	hsa-miR-98	
QM005-02	A04	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM005-02	A05	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM005-02	A06	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM005-02	A07	HmiRQP0408	MIMAT0006764	hsa-miR-320d	
QM005-02	A08	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM005-02	A09	HmiRQP0314	MIMAT0001413	hsa-miR-20b-5p	
QM005-02	A10	HmiRQP0407	MIMAT0005793	hsa-miR-320c	
QM005-02	A11	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM005-02	A12	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM005-02	B01	HmiRQP0159	MIMAT0000691	hsa-miR-130b-3p	
QM005-02	B02	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM005-02	B03	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM005-02	B04	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM005-02	B05	HmiRQP0006	MIMAT0000064	hsa-let-7c	
QM005-02	B06	HmiRQP0004	MIMAT0000063	hsa-let-7b-5p	
QM005-02	B07	HmiRQP0170	MIMAT0000428	hsa-miR-135a-5p	
QM005-02	B08	HmiRQP0255	MIMAT0000072	hsa-miR-18a-5p	
QM005-02	B09	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM005-02	B10	HmiRQP0256	MIMAT0001412	hsa-miR-18b-5p	
QM005-02	B11	HmiRQP0598	MIMAT0002846	hsa-miR-520c-3p	
QM005-02	B12	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM005-02	C01	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM005-02	C02	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM005-02	C03	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM005-02	C04	HmiRQP0286	MIMAT0001080	hsa-miR-196b-5p	
QM005-02	C05	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM005-02	C06	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM005-02	C07	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM005-02	C08	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM005-02	C09	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM005-02	C10	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM005-02	C11	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM005-02	C12	HmiRQP0405	MIMAT0000510	hsa-miR-320a	
QM005-02	D01	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM005-02	D02	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM005-02	D03	HmiRQP0612	MIMAT0005455	hsa-miR-520c-5p	
QM005-02	D04	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM005-02	D05	HmiRQP0299	MIMAT0004571	hsa-miR-200b-5p	
QM005-02	D06	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM005-02	D07	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM005-02	D08	HmiRQP0005	MIMAT0004483	hsa-let-7c*	
QM005-02	D09	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM005-02	D10	HmiRQP0009	MIMAT0004485	hsa-let-7e-3p	
QM005-02	D11	HmiRQP0011	MIMAT0004486	hsa-let-7f-1-3p	
QM005-02	D12	HmiRQP0013	MIMAT0004487	hsa-let-7f-2-3p	
QM005-02	E01	HmiRQP0016	MIMAT0004585	hsa-let-7i-3p	
QM005-02	E02	HmiRQP0019	MIMAT0004512	hsa-miR-100-3p	
QM005-02	E03	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM005-02	E04	HmiRQP0033	MIMAT0004556	hsa-miR-10b-3p	
QM005-02	E05	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM005-02	E06	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM005-02	E07	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM005-02	E08	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM005-02	E09	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM005-02	E10	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM005-02	E11	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM005-02	E12	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM005-02	F01	HmiRQP0304	MIMAT0002810	hsa-miR-202-5p	
QM005-02	F02	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM005-02	F03	HmiRQP0333	MIMAT0004495	hsa-miR-22-5p	
QM005-02	F04	HmiRQP0374	MIMAT0004515	hsa-miR-29b-2-5p	
QM005-02	F05	HmiRQP0376	MIMAT0004673	hsa-miR-29c-5p	
QM005-02	F06	HmiRQP0402	MIMAT0004504	hsa-miR-31-3p	
QM005-02	F07	HmiRQP0422	MIMAT0004703	hsa-miR-335-3p	
QM005-02	F08	HmiRQP0539	MIMAT0004768	hsa-miR-497-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM005-02	F09	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM005-02	F10	HmiRQP1309	MIMAT0010195	hsa-let-7a-2-3p	
QM005-02	F11	HmiRQP1523	MIMAT0015072	hsa-miR-320e	
QM005-02	F12	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM005-02	G01	HmiRQP0097	MIMAT0004603	hsa-miR-125b-2-3p	
QM005-02	G02	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM005-02	G03	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM005-02	G04	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM005-02	G05	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM005-02	G06	HmiRQP0229	MIMAT0000071	hsa-miR-17-3p	
QM005-02	G07	HmiRQP0257	MIMAT0004751	hsa-miR-18b-3p	
QM005-02	G08	HmiRQP0264	MIMAT0001618	hsa-miR-191-3p	
QM005-02	G09	HmiRQP0279	MIMAT0004767	hsa-miR-193b-5p	
QM005-02	G10	HmiRQP0285	MIMAT0004562	hsa-miR-196a-3p	
QM005-02	G11	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM005-02	G12	HmiRQP0294	MIMAT0004491	hsa-miR-19b-1-5p	
QM005-02	H01	NC	NC	NC	
QM005-02	H02	NC	NC	NC	
QM005-02	H03	HmiRQP9001	NR_002752	HK1	
QM005-02	H04	HmiRQP9011	NR_002750	HK2	
QM005-02	H05	HmiRQP9021	NR_002745	HK3	
QM005-02	H06	HmiRQP9051	NR_002746	HK4	
QM005-02	H07	HmiRQP9061	NR_002744	HK5	
QM005-02	H08	HmiRQP9071	NR_002450	HK6	
QM005-02	H09	RT	RT	RT	
QM005-02	H10	RT	RT	RT	
QM005-02	H11	PCR	PCR	PCR	
QM005-02	H12	PCR	PCR	PCR	