

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM050_01	A01	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM050_01	A02	HmiRQP0197	MIMAT0002809	hsa-miR-146b-5p	
QM050_01	A03	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM050_01	A04	HmiRQP0190	MIMAT0000436	hsa-miR-144-3p	
QM050_01	A05	HmiRQP0212	MIMAT0004697	hsa-miR-151a-5p	
QM050_01	A06	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM050_01	A07	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM050_01	A08	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM050_01	A09	HmiRQP0378	MIMAT0000688	hsa-miR-301a-3p	
QM050_01	A10	HmiRQP0193	MIMAT0006789	hsa-miR-1468-5p	
QM050_01	A11	HmiRQP0248	MIMAT0000456	hsa-miR-186-5p	
QM050_01	A12	HmiRQP0540	MIMAT0002820	hsa-miR-497-5p	
QM050_01	B01	HmiRQP0198	MIMAT0004766	hsa-miR-146b-3p	
QM050_01	B02	HmiRQP0317	MIMAT0000267	hsa-miR-210-3p	
QM050_01	B03	HmiRQP0012	MIMAT0000067	hsa-let-7f-5p	
QM050_01	B04	HmiRQP0134	MIMAT0005880	hsa-miR-1290	
QM050_01	B05	HmiRQP0233	MIMAT0004558	hsa-miR-181a-2-3p	
QM050_01	B06	HmiRQP0015	MIMAT0000414	hsa-let-7g-5p	
QM050_01	B07	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM050_01	B08	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM050_01	B09	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM050_01	B10	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM050_01	B11	HmiRQP0399	MIMAT0000693	hsa-miR-30e-3p	
QM050_01	B12	HmiRQP0078	MIMAT0005898	hsa-miR-1246	
QM050_01	C01	HmiRQP0228	MIMAT0004518	hsa-miR-16-2-3p	
QM050_01	C02	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM050_01	C03	HmiRQP1496	MIMAT0015041	hsa-miR-1260b	
QM050_01	C04	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM050_01	C05	HmiRQP2128	MIMAT0019022	hsa-miR-4488	
QM050_01	C06	HmiRQP0344	MIMAT0000078	hsa-mir-23a	
QM050_01	C07	HmiRQP0349	MIMAT0000080	hsa-mir-24-3p	
QM050_01	C08	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM050_01	C09	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM050_01	C10	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM050_01	C11	HmiRQP2147	MIMAT0019041	hsa-miR-4505	
QM050_01	C12	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM050_01	D01	HmiRQP2402	MIMAT0019871	hsa-miR-4741	
QM050_01	D02	HmiRQP0006	MIMAT0000064	hsa-let-7c-5p	
QM050_01	D03	HmiRQP0222	MIMAT0004488	hsa-mir-15a	
QM050_01	D04	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM050_01	D05	HmiRQP0486	MIMAT0003329	hsa-miR-411-5p	
QM050_01	D06	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM050_01	D07	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM050_01	D08	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM050_01	D09	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM050_01	D10	HmiRQP0392	MIMAT0004589	hsa-miR-30b-3p	
QM050_01	D11	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM050_01	D12	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM050_01	E01	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM050_01	E02	HmiRQP1574	MIMAT0016878	hsa-miR-4257	
QM050_01	E03	HmiRQP0394	MIMAT0004674	hsa-miR-30c-1-3p	
QM050_01	E04	HmiRQP0395	MIMAT0004550	hsa-miR-30c-2-3p	
QM050_01	E05	HmiRQP4680	MIMAT0037306	hsa-miR-103a-1-5p	
QM050_01	E06	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM050_01	E07	HmiRQP0023	MIMAT0007402	hsa-miR-103b	
QM050_01	E08	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM050_01	E09	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM050_01	E10	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM050_01	E11	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM050_01	E12	HmiRQP0305	MIMAT0000264	hsa-miR-203a-3p	
QM050_01	F01	HmiRQP2343	MIMAT0019813	hsa-miR-203b-5p	
QM050_01	F02	HmiRQP2344	MIMAT0019814	hsa-miR-203b-3p	
QM050_01	F03	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM050_01	F04	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM050_01	F05	HmiRQP0340	MIMAT0004569	hsa-miR-222-5p	
QM050_01	F06	HmiRQP0313	MIMAT0004752	hsa-mir-20b	
QM050_01	F07	HmiRQP0315	MIMAT0004494	hsa-mir-21	
QM050_01	F08	HmiRQP0276	MIMAT0000459	hsa-mir-193a-3p	
QM050_01	F09	HmiRQP0486	MIMAT0003329	hsa-mir-411	
QM050_01	F10	HmiRQP2626	MIMAT0022259	hsa-mir-5100	
QM050_01	F11	HmiRQP3407	MIMAT0022721	hsa-mir-1247-3p	
QM050_01	F12	HmiRQP0191	MIMAT0004601	hsa-mir-145	
QM050_01	G01	HmiRQP0337	MIMAT0004568	hsa-mir-221	
QM050_01	G02	HmiRQP0339	MIMAT0000279	hsa-mir-222	
QM050_01	G03	HmiRQP0341	MIMAT0004570	hsa-mir-223	
QM050_01	G04	HmiRQP0344	MIMAT0000078	hsa-mir-23a-3p	
QM050_01	G05	HmiRQP0439	MIMAT0004557	hsa-mir-34a	
QM050_01	G06	HmiRQP0504	MIMAT0001541	hsa-mir-449a	
QM050_01	G07	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM050_01	G08	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM050_01	G09	HmiRQP0442	MIMAT0000685	hsa-miR-34b-5p	
QM050_01	G10	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM050_01	G11	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	
QM050_01	G12	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM050_01	H01	NC	NC	NC	
QM050_01	H02	NC	NC	NC	
QM050_01	H03	HKG1	HKG1	HKG1	
QM050_01	H04	HKG2	HKG2	HKG2	
QM050_01	H05	HKG3	HKG3	HKG3	
QM050_01	H06	HKG4	HKG4	HKG4	
QM050_01	H07	HKG5	HKG5	HKG5	
QM050_01	H08	HKG6	HKG6	HKG6	
QM050_01	H09	RT	RT	RT	
QM050_01	H10	RT	RT	RT	
QM050_01	H11	PCR	PCR	PCR	
QM050_01	H12	PCR	PCR	PCR	
QM050_02	A01	HmiRQP0070	MIMAT0005591	hsa-miR-1236	
QM050_02	A02	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM050_02	A03	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM050_02	A04	HmiRQP0097	MIMAT0004603	hsa-miR-125b-2-3p	
QM050_02	A05	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM050_02	A06	HmiRQP1295	MIMAT0009198	hsa-miR-224-3p	
QM050_02	A07	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM050_02	A08	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM050_02	A09	HmiRQP0171	MIMAT0004698	hsa-mir-135b	
QM050_02	A10	HmiRQP0203	MIMAT0004549	hsa-mir-148a	
QM050_02	A11	HmiRQP0358	MIMAT0004501	hsa-mir-27a	
QM050_02	A12	HmiRQP0003	MIMAT0004482	hsa-let-7b	
QM050_02	B01	HmiRQP0254	MIMAT0002891	hsa-mir-18a	
QM050_02	B02	HmiRQP0318	MIMAT0000268	hsa-mir-211	
QM050_02	B03	HmiRQP0318	MIMAT0000268	hsa-mir-211-5p	
QM050_02	B04	HmiRQP2236	MIMAT0019704	hsa-miR-4644	
QM050_02	B05	HmiRQP0031	MIMAT0004555	hsa-mir-10a	
QM050_02	B06	HmiRQP0175	MIMAT0000429	hsa-mir-137	
QM050_02	B07	HmiRQP0211	MIMAT0000757	hsa-mir-151a	
QM050_02	B08	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM050_02	B09	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM050_02	B10	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM050_02	B11	HmiRQP0678	MIMAT0003242	hsa-miR-577	
QM050_02	B12	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM050_02	C01	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM050_02	C02	HmiRQP0297	MIMAT0001620	hsa-miR-200a-5p	
QM050_02	C03	HmiRQP0686	MIMAT0003249	hsa-miR-584-5p	
QM050_02	C04	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM050_02	C05	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM050_02	C06	HmiRQP4536	MIMAT0030019	hsa-miR-7704	
QM050_02	C07	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM050_02	C08	HmiRQP0280	MIMAT0000460	hsa-miR-194-5p	
QM050_02	C09	HmiRQP0322	MIMAT0000272	hsa-miR-215-5p	
QM050_02	C10	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM050_02	C11	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM050_02	C12	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM050_02	D01	HmiRQP0186	MIMAT0000434	hsa-mir-142-3p	
QM050_02	D02	HmiRQP0276	MIMAT0000459	hsa-mir-193a	
QM050_02	D03	HmiRQP2458	MIMAT0019927	hsa-miR-4772-3p	
QM050_02	D04	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM050_02	D05	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM050_02	D06	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM050_02	D07	HmiRQP0630	MIMAT0004806	hsa-miR-548c-5p	
QM050_02	D08	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM050_02	D09	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM050_02	D10	HmiRQP0523	MIMAT0002177	hsa-mir-486-5p	
QM050_02	D11	HmiRQP0832	MIMAT0000092	hsa-mir-92a-3p	
QM050_02	D12	HmiRQP0345	MIMAT0004496	hsa-miR-23a-5p	
QM050_02	E01	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM050_02	E02	HmiRQP1529	MIMAT0015079	hsa-miR-3195	
QM050_02	E03	HmiRQP2192	MIMAT0019337	hsa-miR-3960	
QM050_02	E04	HmiRQP2495	MIMAT0019964	hsa-miR-4792	
QM050_02	E05	HmiRQP3416	MIMAT0022727	hsa-miR-1307-5p	
QM050_02	E06	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	
QM050_02	E07	HmiRQP0809	MIMAT0004949	hsa-miR-877-5p	
QM050_02	E08	HmiRQP0159	MIMAT0000691	hsa-miR-130b-3p	
QM050_02	E09	HmiRQP0491	MIMAT0001340	hsa-miR-423-3p	
QM050_02	E10	HmiRQP2450	MIMAT0019919	hsa-miR-4767	
QM050_02	E11	HmiRQP0835	MIMAT0003218	hsa-miR-92b-3p	
QM050_02	E12	HmiRQP0225	MIMAT0000417	hsa-mir-15b-5p	
QM050_02	F01	HmiRQP0406	MIMAT0005792	hsa-mir-320b	
QM050_02	F02	HmiRQP0172	MIMAT0000758	hsa-miR-135b-5p	
QM050_02	F03	HmiRQP0278	MIMAT0002819	hsa-miR-193b-3p	
QM050_02	F04	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM050_02	F05	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM050_02	F06	HmiRQP0516	MIMAT0003150	hsa-miR-455-5p	
QM050_02	F07	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM050_02	F08	HmiRQP0404	MIMAT0000090	hsa-miR-32-5p	
QM050_02	F09	HmiRQP0786	MIMAT0000252	hsa-miR-7-5p	
QM050_02	F10	HmiRQP0213	MIMAT0000438	hsa-miR-152-3p	
QM050_02	F11	HmiRQP0298	MIMAT0000682	hsa-miR-200a-3p	
QM050_02	F12	HmiRQP0849	MIMAT0004987	hsa-miR-944	
QM050_02	G01	HmiRQP0037	MIMAT0005825	hsa-miR-1180-3p	
QM050_02	G02	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM050_02	G03	HmiRQP0402	MIMAT0004504	hsa-miR-31-3p	
QM050_02	G04	HmiRQP0510	MIMAT0001635	hsa-miR-452-5p	
QM050_02	G05	HmiRQP2779	MIMAT0024599	hsa-mir-6126	
QM050_02	G06	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM050_02	G07	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM050_02	G08	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM050_02	G09	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM050_02	G10	HmiRQP0154	MIMAT0005951	hsa-miR-1307-3p	
QM050_02	G11	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM050_02	G12	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM050_02	H01	NC	NC	NC	
QM050_02	H02	NC	NC	NC	
QM050_02	H03	HKG1	HKG1	HKG1	
QM050_02	H04	HKG2	HKG2	HKG2	
QM050_02	H05	HKG3	HKG3	HKG3	
QM050_02	H06	HKG4	HKG4	HKG4	
QM050_02	H07	HKG5	HKG5	HKG5	
QM050_02	H08	HKG6	HKG6	HKG6	
QM050_02	H09	RT	RT	RT	
QM050_02	H10	RT	RT	RT	
QM050_02	H11	PCR	PCR	PCR	
QM050_02	H12	PCR	PCR	PCR	
QM050_03	A01	HmiRQP4732	MIMAT0037313	hsa-miR-375-5p	
QM050_03	A02	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM050_03	A03	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM050_03	A04	HmiRQP0099	MIMAT0000445	hsa-mir-126-3p	
QM050_03	A05	HmiRQP0210	MIMAT0000451	hsa-mir-150-5p	
QM050_03	A06	HmiRQP0229	MIMAT0000071	hsa-mir-17-3p	
QM050_03	A07	HmiRQP0234	MIMAT0000257	hsa-mir-181b-5p	
QM050_03	A08	HmiRQP0354	MIMAT0000082	hsa-mir-26a-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM050_03	A09	HmiRQP0373	MIMAT0000100	hsa-mir-29b-3p	
QM050_03	A10	HmiRQP2084	MIMAT0018976	hsa-mir-4454	
QM050_03	A11	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM050_03	A12	HmiRQP0517	MIMAT0004761	hsa-miR-483-5p	
QM050_03	B01	HmiRQP0466	MIMAT0000728	hsa-miR-375	
QM050_03	B02	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM050_03	B03	HmiRQP0118	MIMAT0005929	hsa-miR-1275	
QM050_03	B04	HmiRQP0674	MIMAT0004795	hsa-miR-574-5p	
QM050_03	B05	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM050_03	B06	HmiRQP0028	MIMAT0004672	hsa-mir-106b	
QM050_03	B07	HmiRQP0229	MIMAT0000071	hsa-mir-17	
QM050_03	B08	HmiRQP0311	MIMAT0004493	hsa-mir-20a	
QM050_03	B09	HmiRQP0836	MIMAT0004509	hsa-mir-93	
QM050_03	B10	HmiRQP0857	MIMAT0000689	hsa-mir-99b-5p	
QM050_03	B11	HmiRQP0058	MIMAT0005458	hsa-miR-1224-5p	
QM050_03	B12	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM050_03	C01	HmiRQP0183	MIMAT0004598	hsa-mir-141	
QM050_03	C02	HmiRQP0301	MIMAT0004657	hsa-mir-200c	
QM050_03	C03	HmiRQP0434	MIMAT0004692	hsa-miR-340-5p	
QM050_03	C04	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM050_03	C05	HmiRQP0358	MIMAT0004501	hsa-miR-27a-5p	
QM050_03	C06	HmiRQP0264	MIMAT0001618	hsa-miR-191-3p	
QM050_03	C07	HmiRQP0371	MIMAT0000086	hsa-mir-29a-3p	
QM050_03	C08	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM050_03	C09	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM050_03	C10	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM050_03	C11	HmiRQP1314	MIMAT0009197	hsa-miR-205-3p	
QM050_03	C12	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM050_03	D01	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM050_03	D02	HmiRQP0347	MIMAT0004587	hsa-miR-23b-5p	
QM050_03	D03	HmiRQP0325	MIMAT0000274	hsa-miR-217-5p	
QM050_03	D04	HmiRQP0154	MIMAT0005951	hsa-mir-1307	
QM050_03	D05	HmiRQP0161	MIMAT0000426	hsa-mir-132-3p	
QM050_03	D06	HmiRQP0311	MIMAT0004493	hsa-mir-20a-3p	
QM050_03	D07	HmiRQP0364	MIMAT0000690	hsa-mir-296-5p	
QM050_03	D08	HmiRQP0156	MIMAT0000425	hsa-mir-130a	
QM050_03	D09	HmiRQP0546	MIMAT0002872	hsa-mir-501	
QM050_03	D10	HmiRQP0390	MIMAT0000088	hsa-mir-30a-3p	
QM050_03	D11	HmiRQP0446	MIMAT0000703	hsa-mir-361-5p	
QM050_03	D12	HmiRQP0220	MIMAT0004658	hsa-mir-155	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM050_03	E01	HmiRQP0098	MIMAT0000444	hsa-mir-126	
QM050_03	E02	HmiRQP0285	MIMAT0004562	hsa-miR-196a-3p	
QM050_03	E03	HmiRQP0378	MIMAT0000688	hsa-mir-301a	
QM050_03	E04	HmiRQP1554	MIMAT0016858	hsa-miR-4306	
QM050_03	E05	HmiRQP0195	MIMAT0004608	hsa-mir-146a	
QM050_03	E06	HmiRQP0033	MIMAT0004556	hsa-mir-10b	
QM050_03	E07	HmiRQP2171	MIMAT0019064	hsa-miR-4525	
QM050_03	E08	HmiRQP0370	MIMAT0004503	hsa-mir-29a	
QM050_03	E09	HmiRQP0055	MIMAT0004590	hsa-mir-122	
QM050_03	E10	HmiRQP0187	MIMAT0004599	hsa-mir-143	
QM050_03	E11	HmiRQP0782	MIMAT0004284	hsa-miR-675-5p	
QM050_03	E12	HmiRQP1310	MIMAT0006790	hsa-miR-675-3p	
QM050_03	F01	NC	NC	NC	
QM050_03	F02	NC	NC	NC	
QM050_03	F03	HKG1	HKG1	HKG1	
QM050_03	F04	HKG2	HKG2	HKG2	
QM050_03	F05	HKG3	HKG3	HKG3	
QM050_03	F06	HKG4	HKG4	HKG4	
QM050_03	F07	HKG5	HKG5	HKG5	
QM050_03	F08	HKG6	HKG6	HKG6	
QM050_03	F09	RT	RT	RT	
QM050_03	F10	RT	RT	RT	
QM050_03	F11	PCR	PCR	PCR	
QM050_03	F12	PCR	PCR	PCR	