

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM016-01	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1	
QM016-01	A02	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM016-01	A03	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM016-01	A04	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM016-01	A05	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM016-01	A06	HmiRQP0175	MIMAT0000429	hsa-miR-137	
QM016-01	A07	HmiRQP0180	MIMAT0000250	hsa-miR-139-5p	
QM016-01	A08	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM016-01	A09	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM016-01	A10	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM016-01	A11	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM016-01	A12	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM016-01	B01	HmiRQP0459	MIMAT0000724	hsa-miR-372	
QM016-01	B02	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM016-01	B03	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM016-01	B04	HmiRQP0177	MIMAT0000430	hsa-miR-138-5p	
QM016-01	B05	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM016-01	B06	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM016-01	B07	HmiRQP0276	MIMAT0000459	hsa-miR-193a-3p	
QM016-01	B08	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM016-01	B09	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM016-01	B10	HmiRQP0364	MIMAT0000690	hsa-miR-296-5p	
QM016-01	B11	HmiRQP0381	MIMAT0000684	hsa-miR-302a-3p	
QM016-01	B12	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	
QM016-01	C01	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM016-01	C02	HmiRQP0018	MIMAT0000098	hsa-miR-100-5p	
QM016-01	C03	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	
QM016-01	C04	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM016-01	C05	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM016-01	C06	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM016-01	C07	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM016-01	C08	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM016-01	C09	HmiRQP0434	MIMAT0004692	hsa-miR-340-5p	
QM016-01	C10	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM016-01	C11	HmiRQP0305	MIMAT0000264	hsa-miR-203	
QM016-01	C12	HmiRQP0424	MIMAT0004695	hsa-miR-337-5p	
QM016-01	D01	HmiRQP0423	MIMAT0000754	hsa-miR-337-3p	
QM016-01	D02	HmiRQP0125	MIMAT0000424	hsa-miR-128	
QM016-01	D03	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM016-01	D04	HmiRQP0435	MIMAT0004694	hsa-miR-342-5p	
QM016-01	D05	HmiRQP0458	MIMAT0004687	hsa-miR-371a-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM016-01	D06	HmiRQP0452	MIMAT0000719	hsa-miR-367-3p	
QM016-01	D07	HmiRQP0850	MIMAT0000094	hsa-miR-95	
QM016-01	D08	HmiRQP0425	MIMAT0004701	hsa-miR-338-5p	
QM016-01	D09	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM016-01	D10	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM016-01	D11	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM016-01	D12	HmiRQP0426	MIMAT0000763	hsa-miR-338-3p	
QM016-01	E01	HmiRQP0419	MIMAT0004700	hsa-miR-331-5p	
QM016-01	E02	HmiRQP0437	MIMAT0000772	hsa-miR-345-5p	
QM016-01	E03	HmiRQP0457	MIMAT0000723	hsa-miR-371a-3p	
QM016-01	E04	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM016-01	E05	HmiRQP0420	MIMAT0000760	hsa-miR-331-3p	
QM016-01	E06	HmiRQP0418	MIMAT0004693	hsa-miR-330-5p	
QM016-01	E07	HmiRQP0436	MIMAT0000753	hsa-miR-342-3p	
QM016-01	E08	HmiRQP0287	MIMAT0000227	hsa-miR-197-3p	
QM016-01	E09	HmiRQP0365	MIMAT0004679	hsa-miR-296-3p	
QM016-01	E10	HmiRQP0179	MIMAT0004552	hsa-miR-139-3p	
QM016-01	E11	HmiRQP0456	MIMAT0000722	hsa-miR-370	
QM016-01	E12	HmiRQP0417	MIMAT0000751	hsa-miR-330-3p	
QM016-01	F01	HmiRQP0277	MIMAT0004614	hsa-miR-193a-5p	
QM016-01	F02	HmiRQP0463	MIMAT0000727	hsa-miR-374a-5p	
QM016-01	F03	HmiRQP0464	MIMAT0004955	hsa-miR-374b-5p	
QM016-01	F04	HmiRQP0204	MIMAT0000243	hsa-miR-148a-3p	
QM016-01	F05	HmiRQP0383	MIMAT0000715	hsa-miR-302b-3p	
QM016-01	F06	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM016-01	F07	HmiRQP0206	MIMAT0000759	hsa-miR-148b-3p	
QM016-01	F08	HmiRQP0322	MIMAT0000272	hsa-miR-215	
QM016-01	F09	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM016-01	F10	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM016-01	F11	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM016-01	F12	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM016-01	G01	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM016-01	G02	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM016-01	G03	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM016-01	G04	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM016-01	G05	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM016-01	G06	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM016-01	G07	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM016-01	G08	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM016-01	G09	HmiRQP0462	MIMAT0004688	hsa-miR-374a-3p	
QM016-01	G10	HmiRQP0231	MIMAT0000270	hsa-miR-181a-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM016-01	G11	HmiRQP0176	MIMAT0004607	hsa-miR-138-1-3p	
QM016-01	G12	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM016-01	H01	NC	NC	NC	
QM016-01	H02	NC	NC	NC	
QM016-01	H03	HmiRQP9001	NR_002752	HK1	
QM016-01	H04	HmiRQP9011	NR_002750	HK2	
QM016-01	H05	HmiRQP9021	NR_002745	HK3	
QM016-01	H06	HmiRQP9051	NR_002746	HK4	
QM016-01	H07	HmiRQP9061	NR_002744	HK5	
QM016-01	H08	HmiRQP9071	NR_002450	HK6	
QM016-01	H09	RT	RT	RT	
QM016-01	H10	RT	RT	RT	
QM016-01	H11	PCR	PCR	PCR	
QM016-01	H12	PCR	PCR	PCR	