

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM032-01	A01	HmiRQP0044	MIMAT0000416	hsa-miR-1	
QM032-01	A02	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM032-01	A03	HmiRQP0180	MIMAT0000250	hsa-miR-139-5p	
QM032-01	A04	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM032-01	A05	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM032-01	A06	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM032-01	A07	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM032-01	A08	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM032-01	A09	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM032-01	A10	HmiRQP0309	MIMAT0000241	hsa-miR-208a	
QM032-01	A11	HmiRQP0310	MIMAT0004960	hsa-miR-208b	
QM032-01	A12	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM032-01	B01	HmiRQP0349	MIMAT0000080	hsa-miR-24-3p	
QM032-01	B02	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM032-01	B03	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM032-01	B04	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM032-01	B05	HmiRQP0188	MIMAT0000435	hsa-miR-143-3p	
QM032-01	B06	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM032-01	B07	HmiRQP0280	MIMAT0000460	hsa-miR-194-5p	
QM032-01	B08	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM032-01	B09	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM032-01	B10	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM032-01	B11	HmiRQP0543	MIMAT0002870	hsa-miR-499a-5p	
QM032-01	B12	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM032-01	C01	HmiRQP0325	MIMAT0000274	hsa-miR-217	
QM032-01	C02	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM032-01	C03	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM032-01	C04	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM032-01	C05	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM032-01	C06	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM032-01	C07	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM032-01	C08	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM032-01	C09	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM032-01	C10	HmiRQP0212	MIMAT0004697	hsa-miR-151a-5p	
QM032-01	C11	HmiRQP0211	MIMAT0000757	hsa-miR-151a-3p	
QM032-01	C12	HmiRQP0473	MIMAT0000732	hsa-miR-378a-3p	
QM032-01	D01	HmiRQP0542	MIMAT0004772	hsa-miR-499a-3p	
QM032-01	D02	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM032-01	D03	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM032-01	D04	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM032-01	D05	HmiRQP0536	MIMAT0002816	hsa-miR-494	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM032-01	D06	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM032-01	D07	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM032-01	D08	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM032-01	D09	HmiRQP0320	MIMAT0000271	hsa-miR-214-3p	
QM032-01	D10	HmiRQP0522	MIMAT0004762	hsa-miR-486-3p	
QM032-01	D11	HmiRQP0111	MIMAT0000446	hsa-miR-127-3p	
QM032-01	D12	HmiRQP0523	MIMAT0002177	hsa-miR-486-5p	
QM032-01	E01	HmiRQP0179	MIMAT0004552	hsa-miR-139-3p	
QM032-01	E02	HmiRQP0408	MIMAT0006764	hsa-miR-320d	
QM032-01	E03	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM032-01	E04	HmiRQP0407	MIMAT0005793	hsa-miR-320c	
QM032-01	E05	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM032-01	E06	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM032-01	E07	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM032-01	E08	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM032-01	E09	HmiRQP0256	MIMAT0001412	hsa-miR-18b-5p	
QM032-01	E10	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM032-01	E11	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM032-01	E12	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM032-01	F01	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM032-01	F02	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM032-01	F03	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM032-01	F04	HmiRQP0167	MIMAT0000770	hsa-miR-133b	
QM032-01	F05	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM032-01	F06	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM032-01	F07	HmiRQP0405	MIMAT0000510	hsa-miR-320a	
QM032-01	F08	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM032-01	F09	HmiRQP0166	MIMAT0000427	hsa-miR-133a	
QM032-01	F10	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	hsa-miR-199b-3p
QM032-01	F11	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM032-01	F12	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM032-01	G01	HmiRQP1455	MIMAT0014999	hsa-miR-378b	
QM032-01	G02	HmiRQP1523	MIMAT0015072	hsa-miR-320e	
QM032-01	G03	HmiRQP1543	MIMAT0016847	hsa-miR-378c	
QM032-01	G04	HmiRQP2028	MIMAT0010214	hsa-miR-151b	
QM032-01	G05	HmiRQP2029	MIMAT0018926	hsa-miR-378d	
QM032-01	G06	HmiRQP2035	MIMAT0018927	hsa-miR-378e	
QM032-01	G07	HmiRQP2040	MIMAT0018932	hsa-miR-378f	
QM032-01	G08	HmiRQP2046	MIMAT0018937	hsa-miR-378g	
QM032-01	G09	HmiRQP2092	MIMAT0018984	hsa-miR-378h	
QM032-01	G10	HmiRQP2371	MIMAT0019840	hsa-miR-451b	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM032-01	G11	HmiRQP2428	MIMAT0019897	hsa-miR-499b-5p	
QM032-01	G12	HmiRQP2429	MIMAT0019898	hsa-miR-499b-3p	
QM032-01	H01	NC	NC	NC	
QM032-01	H02	NC	NC	NC	
QM032-01	H03	HmiRQP9001	NR_002752	HK1	
QM032-01	H04	HmiRQP9011	NR_002750	HK2	
QM032-01	H05	HmiRQP9021	NR_002745	HK3	
QM032-01	H06	HmiRQP9051	NR_002746	HK4	
QM032-01	H07	HmiRQP9061	NR_002744	HK5	
QM032-01	H08	HmiRQP9071	NR_002450	HK6	
QM032-01	H09	RT	RT	RT	
QM032-01	H10	RT	RT	RT	
QM032-01	H11	PCR	PCR	PCR	
QM032-01	H12	PCR	PCR	PCR	