

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM054_01	A01	HmiRQP0031	MIMAT0004555	hsa-miR-10a-3p	
QM054_01	A02	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM054_01	A03	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM054_01	A04	HmiRQP0678	MIMAT0003242	hsa-miR-577	
QM054_01	A05	HmiRQP0032	MIMAT0000253	hsa-miR-10a-5p	
QM054_01	A06	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM054_01	A07	HmiRQP0297	MIMAT0001620	hsa-miR-200a-5p	
QM054_01	A08	HmiRQP0686	MIMAT0003249	hsa-miR-584-5p	
QM054_01	A09	HmiRQP0134	MIMAT0005880	hsa-miR-1290	
QM054_01	A10	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM054_01	A11	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM054_01	A12	HmiRQP4536	MIMAT0030019	hsa-miR-7704	
QM054_01	B01	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM054_01	B02	HmiRQP0280	MIMAT0000460	hsa-miR-194-5p	
QM054_01	B03	HmiRQP0322	MIMAT0000272	hsa-miR-215-5p	
QM054_01	B04	HmiRQP0239	MIMAT0000259	hsa-miR-182-5p	
QM054_01	B05	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM054_01	B06	HmiRQP0497	MIMAT0001536	hsa-miR-429	
QM054_01	B07	HmiRQP0186	MIMAT0000434	hsa-mir-142-3p	
QM054_01	B08	HmiRQP0276	MIMAT0000459	hsa-mir-193a	
QM054_01	B09	HmiRQP2458	MIMAT0019927	hsa-miR-4772-3p	
QM054_01	B10	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM054_01	B11	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM054_01	B12	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM054_01	C01	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM054_01	C02	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM054_01	C03	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM054_01	C04	HmiRQP0630	MIMAT0004806	hsa-miR-548c-5p	
QM054_01	C05	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM054_01	C06	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM054_01	C07	HmiRQP0221	MIMAT0000646	hsa-mir-155-5p	
QM054_01	C08	HmiRQP0523	MIMAT0002177	hsa-mir-486-5p	
QM054_01	C09	HmiRQP0832	MIMAT0000092	hsa-mir-92a-3p	
QM054_01	C10	HmiRQP0378	MIMAT0000688	hsa-miR-301a-3p	
QM054_01	C11	HmiRQP0345	MIMAT0004496	hsa-miR-23a-5p	
QM054_01	C12	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM054_01	D01	HmiRQP0352	MIMAT0000081	hsa-mir-25-3p	
QM054_01	D02	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM054_01	D03	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM054_01	E01	NC	NC	NC	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM054_01	E02	NC	NC	NC	
QM054_01	E03	HKG1	HKG1	HKG1	
QM054_01	E04	HKG2	HKG2	HKG2	
QM054_01	E05	HKG3	HKG3	HKG3	
QM054_01	E06	HKG4	HKG4	HKG4	
QM054_01	E07	HKG5	HKG5	HKG5	
QM054_01	E08	HKG6	HKG6	HKG6	
QM054_01	E09	RT	RT	RT	
QM054_01	E10	RT	RT	RT	
QM054_01	E11	PCR	PCR	PCR	
QM054_01	E12	PCR	PCR	PCR	