

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM048_01	A01	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM048_01	A02	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	
QM048_01	A03	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM048_01	A04	HmiRQP0494	MIMAT0001341	hsa-miR-424-5p	
QM048_01	A05	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM048_01	A06	HmiRQP0006	MIMAT0000064	hsa-let-7c-5p	
QM048_01	A07	HmiRQP0376	MIMAT0004673	hsa-miR-29c-5p	
QM048_01	A08	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM048_01	A09	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM048_01	A10	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM048_01	A11	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM048_01	A12	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM048_01	B01	HmiRQP0493	MIMAT0004749	hsa-miR-424-3p	
QM048_01	B02	HmiRQP0371	MIMAT0000086	hsa-miR-29a-3p	
QM048_01	B03	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM048_01	B04	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM048_01	B05	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM048_01	B06	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM048_01	B07	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM048_01	B08	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM048_01	B09	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM048_01	B10	HmiRQP0326	MIMAT0004565	hsa-miR-218-1-3p	
QM048_01	B11	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM048_01	B12	HmiRQP0275	MIMAT0004543	hsa-miR-192-3p	
QM048_01	C01	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM048_01	C02	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM048_01	C03	HmiRQP0825	MIMAT0000441	hsa-miR-9-5p	
QM048_01	C04	HmiRQP0517	MIMAT0004761	hsa-miR-483-5p	
QM048_01	C05	HmiRQP0466	MIMAT0000728	hsa-miR-375	
QM048_01	C06	HmiRQP0118	MIMAT0005929	hsa-miR-1275	
QM048_01	C07	HmiRQP0134	MIMAT0005880	hsa-miR-1290	
QM048_01	C08	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM048_01	C09	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM048_01	C10	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM048_01	C11	HmiRQP0300	MIMAT0000318	hsa-miR-200b-3p	
QM048_01	C12	HmiRQP0307	MIMAT0000266	hsa-miR-205-5p	
QM048_01	D01	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	
QM048_01	D02	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM048_01	D03	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM048_01	D04	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM048_01	D05	HmiRQP0315	MIMAT0004494	hsa-miR-21	
QM048_01	D06	HmiRQP2084	MIMAT0018976	hsa-miR-4454	
QM048_01	E01	NC	NC	NC	
QM048_01	E02	NC	NC	NC	
QM048_01	E03	HKG1	HKG1	HKG1	
QM048_01	E04	HKG2	HKG2	HKG2	
QM048_01	E05	HKG3	HKG3	HKG3	
QM048_01	E06	HKG4	HKG4	HKG4	
QM048_01	E07	HKG5	HKG5	HKG5	
QM048_01	E08	HKG6	HKG6	HKG6	
QM048_01	E09	RT	RT	RT	
QM048_01	E10	RT	RT	RT	
QM048_01	E11	PCR	PCR	PCR	
QM048_01	E12	PCR	PCR	PCR	