

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM068-1	A01	HmiRQP0175	MIMAT0000429	hsa-miR-137-3p	
QM068-1	A02	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM068-1	A03	HmiRQP0294	MIMAT0004491	hsa-miR-19b-1-5p	
QM068-1	A04	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM068-1	A05	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM068-1	A06	HmiRQP0187	MIMAT0004599	hsa-miR-143-5p	
QM068-1	A07	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM068-1	A08	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM068-1	A09	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM068-1	A10	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM068-1	A11	HmiRQP0110	MIMAT0004604	hsa-miR-127-5p	
QM068-1	A12	HmiRQP0299	MIMAT0004571	hsa-miR-200b-5p	
QM068-1	B01	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM068-1	B02	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM068-1	B03	HmiRQP0844	MIMAT0004982	hsa-miR-939-5p	
QM068-1	B04	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM068-1	B05	HmiRQP0326	MIMAT0004565	hsa-miR-218-1-3p	
QM068-1	B06	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM068-1	B07	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM068-1	B08	HmiRQP0305	MIMAT0000264	hsa-miR-203a-3p	
QM068-1	B09	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM068-1	B10	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM068-1	B11	HmiRQP0785	MIMAT0004553	hsa-miR-7-1-3p	
QM068-1	B12	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM068-1	C01	HmiRQP0226	MIMAT0004489	hsa-miR-16-1-3p	
QM068-1	C02	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM068-1	C03	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM068-1	C04	HmiRQP0348	MIMAT0000079	hsa-miR-24-1-5p	
QM068-1	C05	HmiRQP0313	MIMAT0004752	hsa-miR-20b-3p	
QM068-1	C06	HmiRQP0030	MIMAT0000104	hsa-miR-107	
QM068-1	C07	HmiRQP0183	MIMAT0004598	hsa-miR-141-5p	
QM068-1	C08	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM068-1	C09	HmiRQP0001	MIMAT0004481	hsa-let-7a-3p	
QM068-1	C10	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM068-1	C11	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM068-1	C12	HmiRQP0536	MIMAT0002816	hsa-miR-494-3p	
QM068-1	D01	HmiRQP0243	MIMAT0004560	hsa-miR-183-3p	
QM068-1	D02	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM068-1	D03	HmiRQP0480	MIMAT0000737	hsa-miR-382-5p	
QM068-1	D04	HmiRQP0782	MIMAT0004284	hsa-miR-675-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM068-1	D05	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM068-1	D06	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM068-1	D07	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM068-1	D08	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM068-1	D09	HmiRQP0423	MIMAT0000754	hsa-miR-337-3p	
QM068-1	D10	HmiRQP0113	MIMAT0005796	hsa-miR-1271-5p	
QM068-1	D11	HmiRQP0546	MIMAT0002872	hsa-miR-501-5p	
QM068-1	D12	HmiRQP0709	MIMAT0003273	hsa-miR-605-5p	
QM068-1	E01	HmiRQP0373	MIMAT0000100	hsa-miR-29b-3p	
QM068-1	E02	HmiRQP2386	MIMAT0019855	hsa-miR-4732-5p	
QM068-1	E03	HmiRQP0460	MIMAT0000725	hsa-miR-373-5p	
QM068-1	E04	HmiRQP0452	MIMAT0000719	hsa-miR-367-3p	
QM068-1	E05	HmiRQP0841	MIMAT0004979	hsa-miR-936	
QM068-1	E06	HmiRQP0242	MIMAT0006767	hsa-miR-1827	
QM068-1	E07	HmiRQP2053	MIMAT0018944	hsa-miR-4429	
QM068-1	E08	HmiRQP2101	MIMAT0018994	hsa-miR-4467	
QM068-1	E09	HmiRQP0329	MIMAT0004567	hsa-miR-219a-1-3p	
QM068-1	E10	HmiRQP0042	MIMAT0005798	hsa-miR-1185-5p	
QM068-1	E11	HmiRQP0169	MIMAT0004595	hsa-miR-135a-3p	
QM068-1	E12	HmiRQP0246	MIMAT0004611	hsa-miR-185-3p	
QM068-1	F01	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM068-1	F02	HmiRQP0309	MIMAT0000241	hsa-miR-208a-3p	
QM068-1	F03	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM068-1	F04	HmiRQP0310	MIMAT0004960	hsa-miR-208b-3p	
QM068-1	F05	HmiRQP0132	MIMAT0005942	hsa-miR-1288-3p	
QM068-1	F06	HmiRQP1523	MIMAT0015072	hsa-miR-320e	
QM068-1	F07	HmiRQP2443	MIMAT0019912	hsa-miR-4763-5p	
QM068-1	F08	HmiRQP2302	MIMAT0019771	hsa-miR-4685-5p	
QM068-1	G01	NC	NC	NC	
QM068-1	G02	NC	NC	NC	
QM068-1	G03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM068-1	G04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM068-1	G05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM068-1	G06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM068-1	G07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM068-1	G08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM068-1	G09	RT	RT	RT	
QM068-1	G10	RT	RT	RT	
QM068-1	G11	PCR	PCR	PCR	
QM068-1	G12	PCR	PCR	PCR	