

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM066-1	A01	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM066-1	A02	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM066-1	A03	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM066-1	A04	HmiRQP0028	MIMAT0004672	hsa-miR-106b-3p	
QM066-1	A05	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM066-1	A06	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM066-1	A07	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM066-1	A08	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM066-1	A09	HmiRQP0003	MIMAT0004482	hsa-let-7b-3p	
QM066-1	A10	HmiRQP0280	MIMAT0000460	hsa-miR-194-5p	
QM066-1	A11	HmiRQP0358	MIMAT0004501	hsa-miR-27a-5p	
QM066-1	A12	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM066-1	B01	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM066-1	B02	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM066-1	B03	HmiRQP0170	MIMAT0000428	hsa-miR-135a-5p	
QM066-1	B04	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM066-1	B05	HmiRQP3046	MIMAT0000250	hsa-miR-139-5p	
QM066-1	B06	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM066-1	B07	HmiRQP0317	MIMAT0000267	hsa-miR-210-3p	
QM066-1	B08	HmiRQP0222	MIMAT0004488	hsa-miR-15a-3p	
QM066-1	B09	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM066-1	B10	HmiRQP0177	MIMAT0000430	hsa-miR-138-5p	
QM066-1	B11	HmiRQP0235	MIMAT0000258	hsa-miR-181c-5p	
QM066-1	B12	HmiRQP0006	MIMAT0000064	hsa-let-7c-5p	
QM066-1	C01	HmiRQP0764	MIMAT0004814	hsa-miR-654-3p	
QM066-1	C02	HmiRQP0245	MIMAT0000454	hsa-miR-184	
QM066-1	C03	HmiRQP0047	MIMAT0005865	hsa-miR-1202	
QM066-1	C04	HmiRQP0059	MIMAT0005572	hsa-miR-1225-5p	
QM066-1	C05	HmiRQP0738	MIMAT0003298	hsa-miR-629-3p	
QM066-1	C06	HmiRQP0040	MIMAT0005828	hsa-miR-1183	
QM066-1	C07	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM066-1	C08	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM066-1	C09	HmiRQP0209	MIMAT0004610	hsa-miR-150-3p	
QM066-1	C10	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM066-1	C11	HmiRQP0435	MIMAT0004694	hsa-miR-342-5p	
QM066-1	C12	HmiRQP0522	MIMAT0004762	hsa-miR-486-3p	
QM066-1	D01	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM066-1	D02	HmiRQP0375	MIMAT0000681	hsa-miR-29c-3p	
QM066-1	D03	HmiRQP0203	MIMAT0004549	hsa-miR-148a-5p	
QM066-1	D04	HmiRQP0118	MIMAT0005929	hsa-miR-1275	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM066-1	D05	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM066-1	D06	HmiRQP0401	MIMAT0000089	hsa-miR-31-5p	
QM066-1	D07	HmiRQP0421	MIMAT0000765	hsa-miR-335-5p	
QM066-1	D08	HmiRQP0686	MIMAT0003249	hsa-miR-584-5p	
QM066-1	D09	HmiRQP0214	MIMAT0000439	hsa-miR-153-3p	
QM066-1	D10	HmiRQP0620	MIMAT0003340	hsa-miR-542-5p	
QM066-1	D11	HmiRQP0438	MIMAT0000773	hsa-miR-346	
QM066-1	D12	HmiRQP0526	MIMAT0002804	hsa-miR-488-5p	
QM066-1	E01	HmiRQP0519	MIMAT0002174	hsa-miR-484	
QM066-1	E02	HmiRQP0318	MIMAT0000268	hsa-miR-211-5p	
QM066-1	E03	HmiRQP0717	MIMAT0003281	hsa-miR-613	
QM066-1	E04	HmiRQP0238	MIMAT0000260	hsa-miR-182-3p	
QM066-1	E05	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM066-1	E06	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM066-1	E07	HmiRQP0073	MIMAT0004591	hsa-miR-124-5p	
QM066-1	E08	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM066-1	E09	HmiRQP0370	MIMAT0004503	hsa-miR-29a-5p	
QM066-1	E10	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM066-1	E11	HmiRQP0719	MIMAT0004804	hsa-miR-615-5p	
QM066-1	E12	HmiRQP0459	MIMAT0000724	hsa-miR-372-3p	
QM066-1	F01	HmiRQP0125	MIMAT0000424	hsa-miR-128-3p	
QM066-1	F02	HmiRQP0173	MIMAT0000448	hsa-miR-136-5p	
QM066-1	G01	NC	NC	NC	
QM066-1	G02	NC	NC	NC	
QM066-1	G03	hsnRNA U6	NR_002752	HK1(RNU6-2)	
QM066-1	G04	hsnoRNA U44	NR_002750	HK2(SNORD44)	
QM066-1	G05	hsnoRNA U48	NR_002745	HK3(SNORD48)	
QM066-1	G06	HsnoRNA U47	NR_002746	HK4(SNORD47)	
QM066-1	G07	HsnoRNA U49	NR_002744	HK5(SNORD49A)	
QM066-1	G08	HsnRNA U68	NR_002450	HK6(SNORD68)	
QM066-1	G09	RT	RT	RT	
QM066-1	G10	RT	RT	RT	
QM066-1	G11	PCR	PCR	PCR	
QM066-1	G12	PCR	PCR	PCR	