

miR-9-1	miR-10b	miR-17-3p	miR-22	miR-23a	miR-24	miR-26a	miR-26b	miR-27a	miR-27b
miR-29a	miR-29b	miR-29c	miR-30a-3p	miR-30a-5p	miR-30b	miR-30c	miR-92	miR-92b	miR-93
miR-95	miR-101-1	miR-103	miR-106a	miR-106b	miR-107	miR-128a	miR-128b	miR-132	miR-134
miR-135b	miR-136	miR-137	miR-140	miR-141	miR-142-3p	miR-149	miR-150	miR-151	miR-153
miR-154	miR-181d	miR-183	miR-185	miR-186	miR-188	miR-190	miR-191	miR-196a	miR-196b
miR-197	miR-198	miR-200b	miR-202	miR-203	miR-205	miR-210	miR-214	miR-215	miR-218
miR-219	miR-221	miR-222	miR-296	miR-372	miR-373	miR-488	miR-100	miR-127	miR-142-5p
miR-31	miR-213	RNU48							

Note: Each miRNA is duplicate spotted