

Let-7a	Let-7b	Let-7c	Let-7d	Let-7e	Let-7f	Let-7g	Let-7i	miR-1	miR-7
miR-9	miR-10a	miR-15a	miR-15b	miR-16	miR-17-5p	miR-18a	miR-18b	miR-19a	miR-19b
miR-20a	miR-21	miR-25	miR-28	miR-34a	miR-99a	miR-122a		miR-125a	miR-125b
miR-126		miR-133a	miR-133b	miR-143	miR-145	miR-146a	miR-146b	miR-148a	miR-155
miR-181a	miR-181b	miR-181c	miR-182	miR-192	miR-194	miR-195	miR-199a		
miR-200a	miR-200c	miR-204	miR-206	miR-216	miR-223	miR-224	miR-342		miR-375
miR-9-1	miR-10b		miR-22	miR-23a	miR-24	miR-26a	miR-26b	miR-27a	miR-27b
miR-29a	miR-29b	miR-29c		miR-30a-5p	miR-30b	miR-30c	miR-92	miR-92b	miR-93
	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-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-200b	miR-202	miR-203	miR-205	miR-210	miR-214	miR-215	miR-218
miR-219	miR-221	miR-222	miR-296				miR-100	miR-127	miR-142-5p
miR-31	miR-213	U6							

Note: Each miRNA is duplicate spotted