

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-124a	miR-125a	miR-125b
miR-126	miR-131	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-199b	miR-199a*
miR-200a	miR-200c	miR-204	miR-206	miR-216	miR-223	miR-224	miR-342	miR-368	miR-375

Note: Each miRNA is duplicate spotted