

Let-7a	Let-7b	Let-7c	Let-7d	Let-7e	Let-7f	Let-7g	Let-7i	miR-1	miR-7
miR-10a	miR-15a	miR-15b	miR-16	miR-21	miR-25	miR-28	miR-133b	miR-145	miR-148
miR-155	miR-182	miR-204	miR-216	miR-224	miR-368	miR-24	miR-101	miR-142	miR-150
miR-151	miR-153	miR-188	miR-196a	miR-197	miR-202	miR-210	miR-214	miR-218	miR-296
miR-144	miR-152	miR-184	miR-337	miR-338	miR-345	miR-371	miR-96	miR-369-5p	miR-369-3p
miR-193a	miR-193b	RUN48							

Note: Each miRNA is duplicate spotted