






## ***FlexmiR™ Analysis using Bio-Plex® Manager™ Software***

FlexmiR MicroRNA analysis kits are compatible with Bio-Plex Instruments and Bio-Plex Manager Software. Calibrate instrument using the standard procedures for Bio-Plex calibration.

### **Human microRNA Analysis**

For analysis of FlexmiR Human Pools 1-5 use the following protocol (.pbx) files.








-  FlexmiR MicroRNA Human Panel-Pool 1.pbx
-  FlexmiR MicroRNA Human Panel-Pool 2.pbx
-  FlexmiR MicroRNA Human Panel-Pool 3.pbx
-  FlexmiR MicroRNA Human Panel-Pool 4.pbx
-  FlexmiR MicroRNA Human Panel-Pool 5.pbx

Refer to FlexmiR Human miRNA Sequence spreadsheet for sequence information.

### **Mouse or Rat microRNA analysis**

For evaluation of Mouse or Rat RNA you should purchase the 2 additional FlexmiR Mouse/Rat Pools. All seven pools are required for the most comprehensive coverage.

For analysis use the following protocol files.

- |  |                    |
|--|--------------------|
|  FlexmiR MicroRNA Human Panel-Pool 1.pbx      | “Universal” Pool 1 |
|  FlexmiR MicroRNA Human Panel-Pool 2.pbx      | “Universal” Pool 2 |
|  FlexmiR MicroRNA Human Panel-Pool 3.pbx      | “Universal” Pool 3 |
|  FlexmiR MicroRNA Human Panel-Pool 4.pbx      | “Universal” Pool 4 |
|  FlexmiR MicroRNA Human Panel-Pool 5.pbx      | “Universal” Pool 5 |
|  |                    |
|  FlexmiR MicroRNA Rodent Panel-Pool 1 (6).pbx | Or Pool 6          |
|  FlexmiR MicroRNA Rodent Panel-Pool 2 (7).pbx | Or Pool 7          |

Refer to FlexmiR Mouse-Rat miRNA Sequence spreadsheet for sequence information.

Mouse Rat Pool1 Referred to as Pool 6 in some Luminex Documentation and IS Templates  
Mouse Rat Pool2 Referred to as Pool 7 in some Luminex Documentation and IS Templates