

Product datasheet

FirePlex[®] miRNA Assay Core Reagent Kit V2 - Biofluid ab218342

FirePlex[®]

2 References

Overview

Product name	FirePlex [®] miRNA Assay Core Reagent Kit V2 - Biofluid
Assay type	Enzyme activity
Assay time	4h 00m
Product overview	FirePlex [®] miRNA Assay Core Reagent Kit V2 - Biofluid (ab218342) contains qualified reagents for the use of our FirePlex [®] miRNA multiplex assays which detect up to 65 literature curated miRNA in only 10 µL of biofluid samples or 100 pg purified total RNA, without the need for RNA purification.

FirePlex[®] miRNA Assay Core Reagent Kit V2 - Biofluid (ab218342) requires the purchase of either:

- Oncology miRNA Panel ([ab218367](#))
- Cardiology miRNA Panel ([ab218368](#))
- Immunology miRNA Panel ([ab218369](#))
- Liver Tox miRNA Panel ([ab218370](#))
- Neurology miRNA Panel ([ab218371](#))
- Kidney Toxicity miRNA Panel ([ab219508](#))
- Cardiac Toxicity miRNA Panel ([ab219509](#))
- Pancreatic Toxicity miRNA Panel ([ab219510](#))

NOTE: FirePlex miRNA panels require the purchase of a Run Buffer. Use our [FirePlex Cytometer Setup Kit \(ab211043\)](#) to select the most appropriate Run Buffer for your flow cytometer.

The assay provides PCR sensitivity while eliminating the need for separate reverse transcription reactions and mitigating amplification biases introduced by target-specific qPCR. This is made possible by combining uniquely encoded FirePlex[®] hydrogel particles with single-step RT-PCR signal amplification using universal primers. The entire assay is performed in 96-well plate format such that users detect up to 65 microRNA targets for each of 96 samples in less than a day's work.

The FirePlex[®] miRNA Assay reliably detects as few as 1000 microRNA copies per sample with a

linear dynamic range of ~5 logs. In addition to increased sensitivity, the assay's multiplexed detection conserves precious sample by detecting multiple microRNA targets from only 10 μ L of biofluid, and has been validated for numerous sample types including plasma, serum, exosomes, urine and saliva.

For a full list of validated sample types, and to learn more about the performance and requirements of these assays, visit <https://www.abcam.com/nav/multiplex-mirna-assays>.

Our integrated free FirePlex Analysis Workbench software enables easy and rapid data analysis, visualization, and export in under ten minutes, and includes key features such as standard curve analysis and publication-quality heat maps and graphs.

Notes	FirePlex [®] is a registered trade mark in the United States and is an unregistered trade mark elsewhere.
Platform	Multiplex

Properties

Storage instructions Please refer to protocols.

Components	1 x 96 tests
10x Rinse Buffer A	1 x 30ml
1x Hyb Buffer	1 x 25ml
2x Rinse Buffer B	1 x 33ml
50x Labeling Buffer	1 x 200 μ l
5x Reporter Solution	1 x 4ml
Catch Plate	1 unit
Digest Buffer	1 x 4ml
dNTP mix	1 x 120 μ l
Filter Plate	1 unit
FirePlex Plate Seal	1 unit
Labeling Diluent	1 x 30ml
Labeling Enzyme	1 x 50 μ l
PCR Buffer	1 x 2ml
PCR Enzyme	1 x 60 μ l
Primer Mix	1 x 250 μ l

Components	1 x 96 tests
Protease Mix	1 x 500µl
RNA Control	1 x 50µl

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors