

Product datasheet

Fructose Assay Kit (Fluorometric) ab241022

2 Images

Overview

<b>Product name</b>	Fructose Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Product overview</b>	Fructose Assay Kit (Fluorometric) (ab241022) measures free fructose that is enzymatically processed with the formation of a metabolite which reacts with the Probe to generate fluorescence (Ex/Em = 535/587 nm). The kit provides a simple, highly sensitive, reliable method suitable for high throughput assay of D-fructose. Glucose interference can be removed by using the Sample Cleanup Mix. The kit can detect fructose in the range of 5 to 500 picomoles/well.

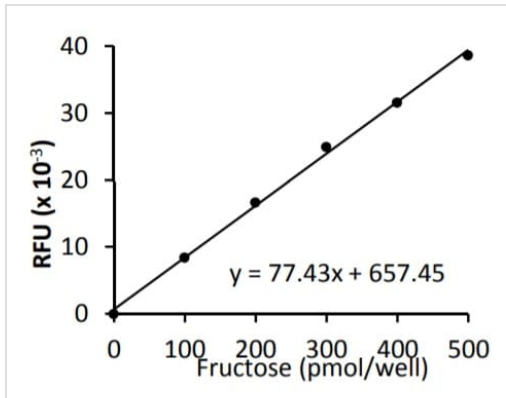
Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	100 tests
Conversion Enzyme	1 vial
Fructose Assay Buffer	1 x 25ml
Fructose Enzyme Mix	1 vial
Fructose Standard (100 mM)	1 x 100µl
Fructose Substrate Mix	1 vial
Probe	1 x 200µl
Sample Cleanup Mix	1 vial

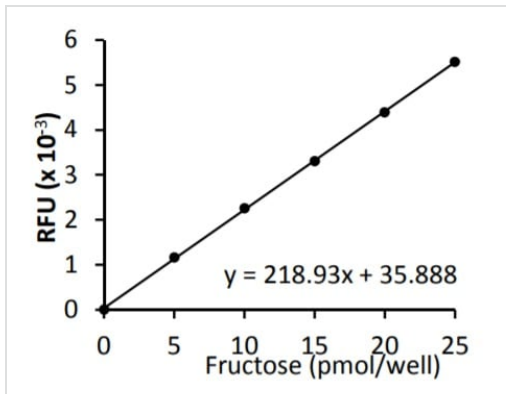
**Relevance** Fructose is a monosaccharide found in many foods and is one of the three most important blood sugars along with glucose and galactose. Fructose is the sweetest naturally occurring sugar, estimated to be twice as sweet as sucrose.

Images



Fructose standard curve generated using this kit protocol.

Fructose standard curve.



Fructose standard curve generated using this kit protocol.

Fructose standard curve.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### 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