

Fructose Assay Kit ab83380

10 References **画像数 4**

製品の概要

製品名	Fructose Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell culture extracts, Tissue Extracts
アッセイタイプ	Quantitative
全工程の試験時間	1h 00m
製品の概要	In Abcam's Fructose Assay Kit, free fructose is enzymatically converted to β-glucose, which is then specifically converted to a product that reacts with OxiRed Probe to generate color (λ=570nm) and fluorescence (Ex/Em=535/587nm). The kit provides a rapid, simple, sensitive, and reliable method suitable for high throughput assay of D-fructose. Do not use this kit with urine, serum, or plasma samples.

Visit our [FAQs page](#) for tips and troubleshooting.

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K619 Fructose Colorimetric/Fluorometric Assay Kit. K619-100 is the same size as the 100 test size of ab83380.</p> <p>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.</p>
------	---

試験プラットフォーム	Microplate reader
------------	-------------------

製品の特性

保存方法	Store at -20°C. Please refer to protocols.
------	--

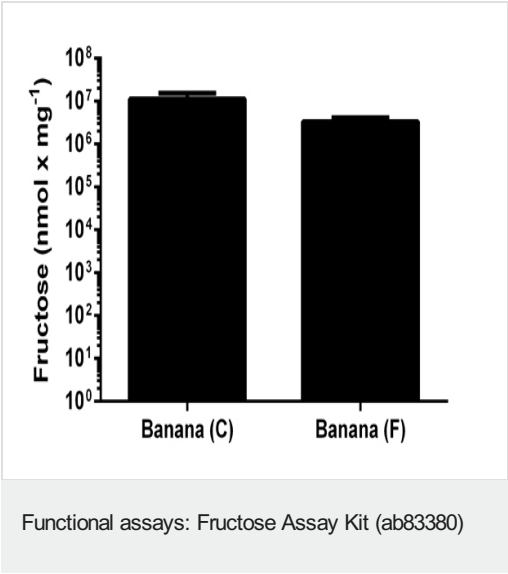
内容	100 tests
Development Enzyme Mix II	1 vial
Fructose Assay Buffer	1 x 25ml
Fructose Converter Mix	1 vial

内容	100 tests
Fructose Standard	1 x 100µl
OxiRed Probe	1 x 200µl

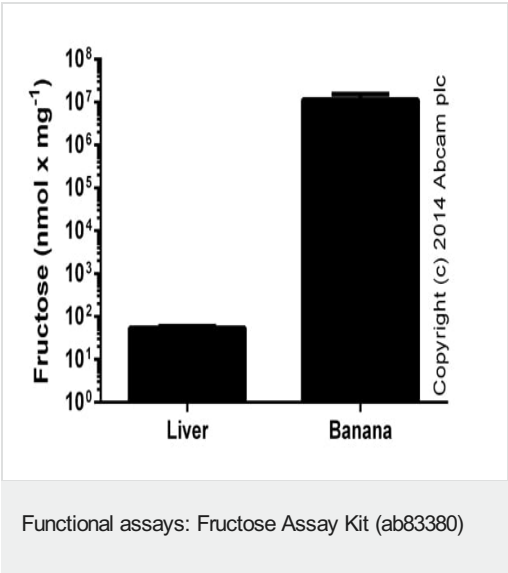
関連性

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.

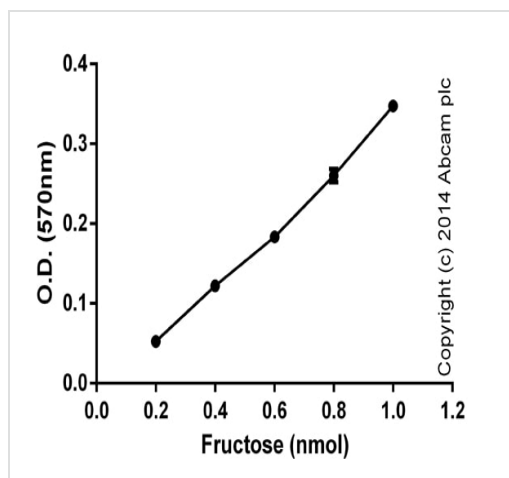
画像



Fructose levels banana, determined using the colorimetric (C) and the fluorometric (F) method (presented per mg of protein) (duplicates; +/- SD).

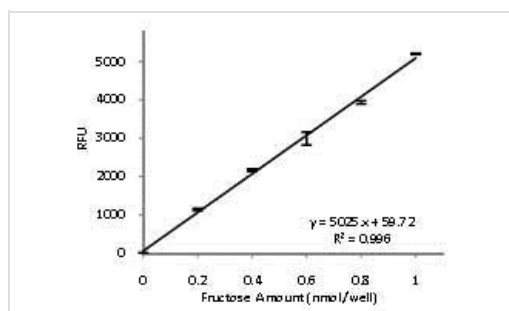


Fructose levels in mouse liver and banana, determined using the colorimetric method and presented per mg of protein (duplicates; +/- SD).



Titration of the fructose (background subtracted, duplicates; +/- SD).

Functional assays: Fructose Assay Kit (ab83380)



Fluorometric Fructose Standard Curve performed following the ab83380 kit protocol.

Functional Studies - Fructose Assay Kit (ab83380)

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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