

# Galactose Assay Kit ab83382

**5 References**   [画像数 5](#)

### 製品の概要

製品名	Galactose Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
検出感度	< 10 nmol/well
全工程の試験時間	0h 40m
製品の概要	Galactose Assay Kit (Colorimetric/Fluorometric) (ab83382) provides a simple and convenient method for direct measurement of galactose levels in various biological samples such as serum, plasma and other body fluids, as well as food and growth media. In this assay, oxidation of galactose generates a product that reacts with the galactose probe to produce color (OD570nm) or fluorescence (Ex/Em 535/587 nm). Liquid samples can be tested directly without purification. The assay is fast, convenient and sensitive. It can also be used as a high throughput assay.

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 K621 Galactose Colorimetric/Fluorometric Assay Kit. K621-100 is the same size as the 100 test size of ab83382.</p> <p>Galactose (C<sub>6</sub>H<sub>12</sub>O<sub>6</sub> MW: 180.16 g/mol) is one of the major naturally occurring sugars, a monosaccharide that differs from glucose only by the configuration of the hydroxyl group at the carbon-4 position. Often present as an anomeric mixture of α-D-galactose and β-D-galactose, this monosaccharide exists abundantly in milk, dairy products and many other food types such as fruits and vegetables.</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.
------	--

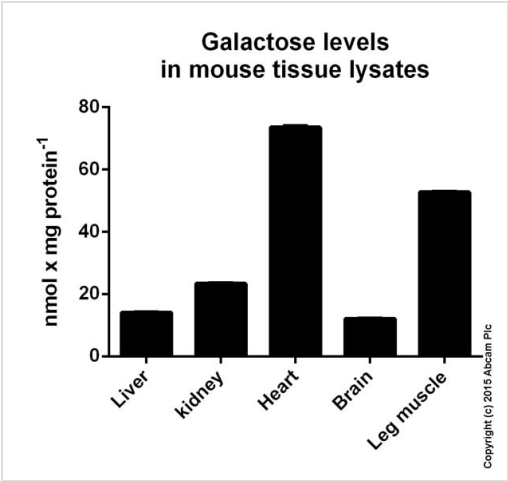
内容	100 tests
Assay Buffer II	1 x 25ml

内容	100 tests
Developer Solution V	1 vial
Galactose Enzyme Mix	1 vial
Galactose Standard	1 x 100µl
OxiRed Probe	1 x 0.2ml

関連性

Galactose (C<sub>6</sub>H<sub>12</sub>O<sub>6</sub> FW: 180.16) is one the major naturally occurring sugars, involved in many biological processes.

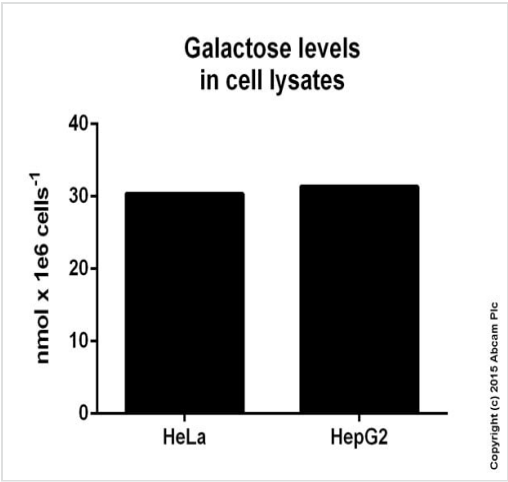
画像



Functional Studies - Galactose Assay Kit (ab83382)

Galactose measured in tissue lysates showing quantity (nmol) per mg of extracted protein.

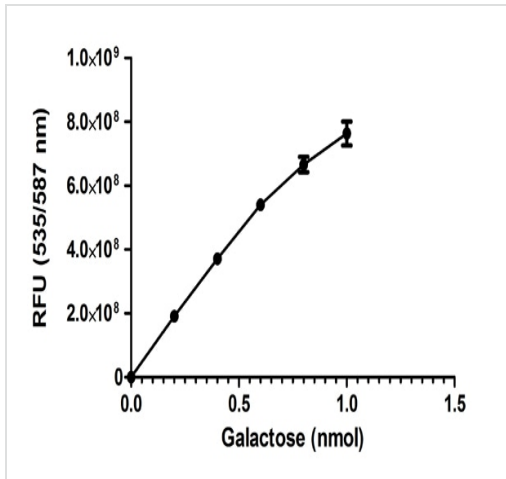
Protein concentration for samples varied from 7 mg/mL to 17 mg/mL. Samples were tested undiluted.



Functional Studies - Galactose Assay Kit (ab83382)

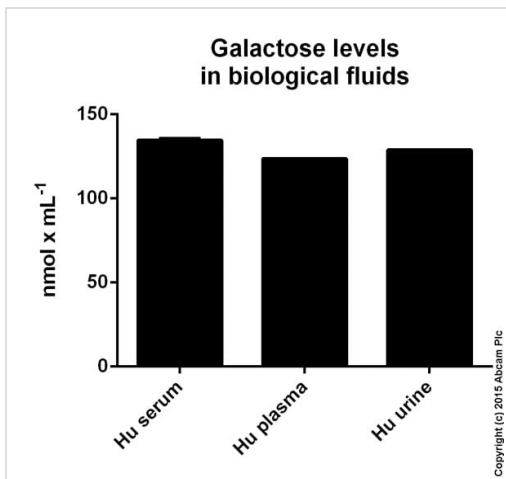
Galactose measured in cell lysates showing quantity (nmol) per 1 mln cells.

Samples with the concentration of 4e6 cells/mL were used. Samples were tested undiluted.



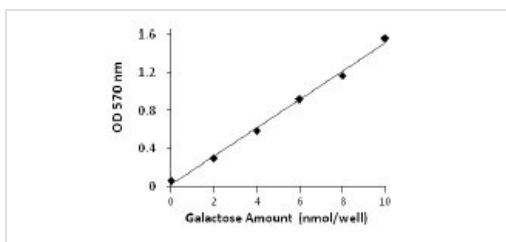
Fluorometric standard curve: mean of duplicates (+/- SD) with background reads subtracted.

Functional Studies - Galactose Assay Kit (ab83382)



Galactose levels measured in biological fluids showing quantity (nmol) per mL. Samples tested undiluted.

Functional Studies - Galactose Assay Kit (ab83382)



Example of Galactose standard curve obtained with ab83382

Functional Studies - Galactose Assay Kit (ab83382)

**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