

Lactose Assay Kit ab83384

5 References 画像数 1

製品の概要

製品名	Lactose Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Milk, Other biological fluids
アッセイタイプ	Quantitative
全工程の試験時間	1h 00m
製品の概要	<p>In Abcam's Lactose Assay Kit Lactose is hydrolyzed to glucose and galactose. The galactose is subsequently oxidized generating color (OD570nm) and fluorescence (Ex/Em 535/587 nm). Free galactose can be corrected by a background control in the absence of lactase. The Lactose Assay Kit provides a simple, convenient, and sensitive means for direct measurement of lactose levels in various biological samples (serum, plasma, other body fluids, food, growth media, etc.). Pretreatment of samples is not required. The kit can be used as a high throughput assay. Visit our FAQs page for tips and troubleshooting.</p>
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K624 Lactose Colorimetric/Fluorometric Assay Kit. K624-100 is the same size as the 100 test size of ab83384.</p> <p>Lactose (C₁₂H₂₂O₁₁ FW: 342.3) is an important naturally occurred disaccharide, consisting of one galactose and one glucose. Milk contains ~2-8% lactose. Some people, particularly infants, lack the enzyme necessary to digest galactose leading to galactose accumulation in blood (Galactosemia) causing enlarged liver, renal failure, cataracts and brain damage.</p>
試験プラットフォーム	Microplate reader

製品の特性

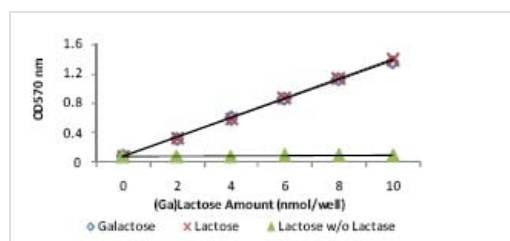
保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
Assay Buffer II	1 x 25ml
Developer Solution V	1 vial
Lactase Enzyme	1 vial

内容	100 tests
Galactose Enzyme Mix	1 vial
Lactose Standard	1 x 100µl
OxiRed Probe	1 x 200µl

関連性 Lactose (C₁₂H₂₂O₁₁ FW: 342.3) is an important naturally occurring disaccharide along with sucrose and maltose. Lactose makes up around 2-8% of milk. It consists of one galactose and one glucose. Some people, particularly infants, lack the enzyme necessary to digest galactose leading to galactose accumulation in blood (Galactosemia) causing enlarged liver, renal failure, cataracts and brain damage.

画像



Lactose Standard Curve: Assays were performed following the kit instructions. The kit detected galactose and lactose equally. In the absence of Lactase, the kit detected galactose, but not lactose.

Functional Studies - Lactose Assay Kit (ab83384)

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