

Lactose Assay Kit ab83384

[5 References](#) [1 图像](#)

概述

产品名称	Lactose Assay试剂盒
检测方法	Colorimetric/Fluorometric
样品类型	Milk, Other biological fluids
检测类型	Quantitative
检测时间	1h 00m
产品概述	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.

说明

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.

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.

平台 Microplate reader

性能

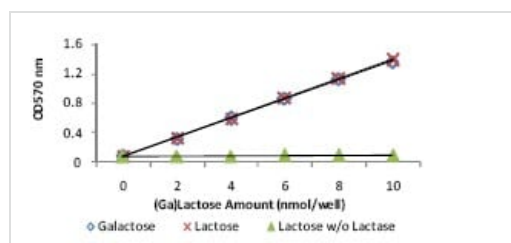
存放说明 Store at -20°C. Please refer to protocols.

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

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

相关性 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.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

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