

Product datasheet

Intestinal Permeability Assay Kit ab272540

2 Images

Overview

Product name	Intestinal Permeability Assay Kit
Detection method	Colorimetric
Sample type	Urine
Assay type	Quantitative
Product overview	<p>Intestinal Permeability Assay Kit (ab272540) is based on measuring the ratio of the absorption of two non-metabolized sugars through the intestines. Test subjects drink a prescribed amount of lactulose and mannitol and the % absorption of these sugars is determined by the amount of excreted lactulose and mannitol measured during the first 6 hours after ingestion. The degree of intestinal permeability is reflected by the ratio of the % absorption of lactulose to % absorption of mannitol. An increase in this ratio indicates increased intestinal permeability since lactulose is only absorbed through intercellular spaces. Lactulose and mannitol are measured in separate assays using the included Lactulose Assay Kit and Mannitol Assay Kit, respectively.</p>

Simple and convenient: Both assays require addition of single working reagent and can be completed within 60 minutes. Both assays are performed at room temperature. No 37°C heater is needed.

High-throughput: Can be readily automated as a high-throughput 96- well plate assay for thousands of samples per day.

Detection Limit: Mannitol - 7 µM; Lactulose - 3 µM.

Notes

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Tested applications **Suitable for:** Functional Studies

Platform Microplate

Properties

Storage instructions

Store at -20°C. Please refer to protocols.

Components	100 tests
Lactulose Assay Buffer	1 x 6ml
Lactulose Enzyme A	1 vial
Lactulose Enzyme B	1 x 120µl
Lactulose Enzyme Buffer	1 x 150µl
Lactulose PMS Solution	1 x 1.5ml
Lactulose Standard	1 x 400µl
Mannitol Standard	1 x 500µl
Mannitol Assay Buffer	1 x 10ml
Mannitol Enzyme	1 x 120µl

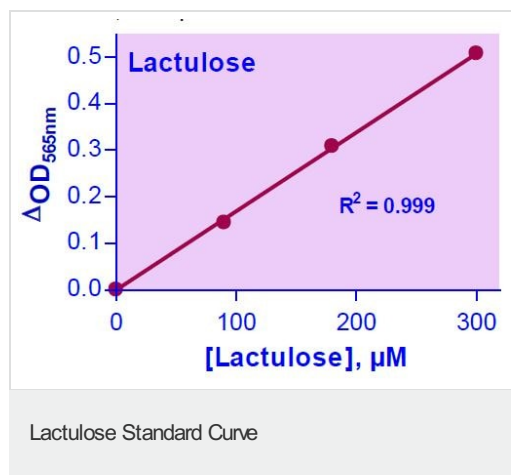
Applications

Our [Abpromise guarantee](#) covers the use of **ab272540** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

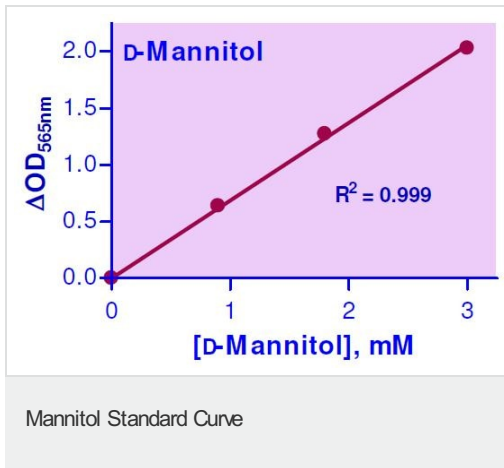
Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

Images



Lactulose Standard Curve

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

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