

alpha-L-Fucosidase Assay Kit ab272525

画像数 2

製品の概要

製品名	alpha-L-Fucosidase Assay Kit
検出方法	Colorimetric
サンプルの種類	Serum, Plasma, Cell Lysate, Tissue Lysate
アッセイタイプ	Quantitative
検出範囲	1 mU/ml - 100 mU/ml
製品の概要	<p>alpha-L-Fucosidase Assay Kit (ab272525) is a non-radioactive, colorimetric assay based on the cleavage of 4-nitrophenol from the synthetic substrate. Nitrophenol becomes intensely colored after addition of the stop reagent. The increase in absorbance at 405 nm after addition of the stop reagent is directly proportional to the enzyme activity.</p> <p>High sensitivity and wide linear range. Linear detection range (10 µL sample): 1 to 100 U/L for a 20 minute reaction.</p> <p>Homogeneous and simple procedure. Simple add-mix-read procedure allows reliable quantitation of fucosidase activity within 20 minutes.</p> <p>Robust and amenable to HTS. All reagents are compatible with highthroughput liquid handling instruments. Can be readily automated to measure thousands of samples per day.</p>
アプリケーション	適用あり: Functional Studies
試験プラットフォーム	Microplate

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

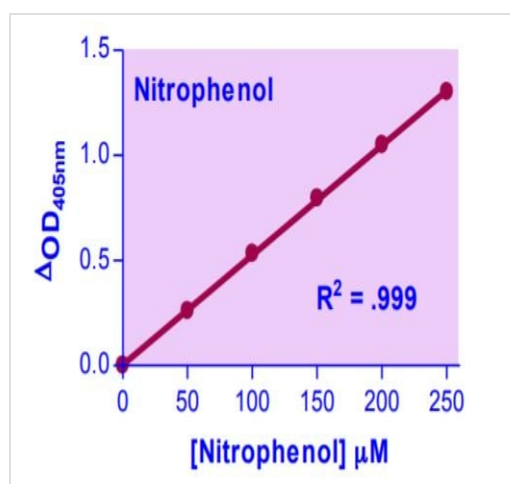
内容	100 tests
Standard (12.5 mM Nitrophenol)	1 x 1ml
Stop Reagent	1 x 12ml
Substrate Buffer	1 x 10ml

アプリケーション

The Abpromise guarantee Abpromise保証は、次のテスト済みアプリケーションにおけるab272525の使用に適用されます。アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

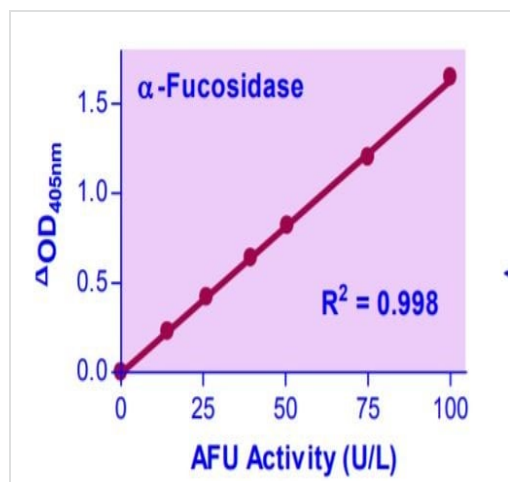
画像



Standard Curve

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

Example Standard Curve.



Activity Titration Curve

Example activity titration curve.

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