

Lipase Assay Kit II (Colorimetric) ab102525

[1 References](#) [2 Images](#)

Overview

Product name	Lipase Assay Kit II (Colorimetric)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
Assay type	Enzyme activity
Assay time	1h 30m
Product overview	Lipase Assay Kit II (Colorimetric) (ab102525) provides a simple, sensitive, and reliable assay for rapid analysis of Lipase activity.

In the lipase assay, Lipases hydrolyze a specific substrate to generate a product which reacts with the DTNB probe to generate color (Absorbance=412 nm). This kit is suitable for high throughput analysis (HTP).

Lipase assay protocol summary:

- add samples and standards to wells
- add reaction mix
- analyze with a microplate reader at 3 min and at between 60 and 90 min

Notes This product is manufactured by BioVision, an Abcam company and was previously called K723 Lipase Activity Colorimetric Assay Kit II. K723-100 is the same size as the 100 test size of ab102525.

Platform Microplate reader

Properties

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

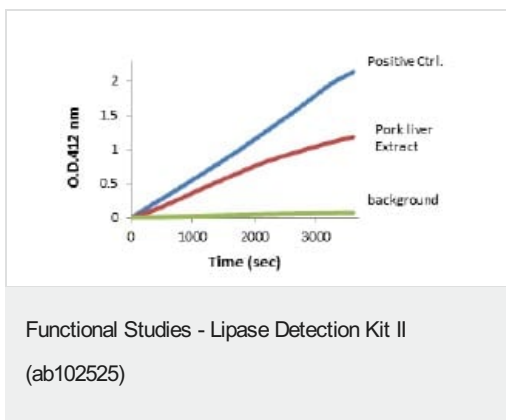
Components	100 tests
Assay Buffer V	1 x 25ml
DTNB	1 vial
Lipase Positive Control	1 vial

Components	100 tests
Lipase Substrate II	1 x 500µl
TNB Standard	1 vial

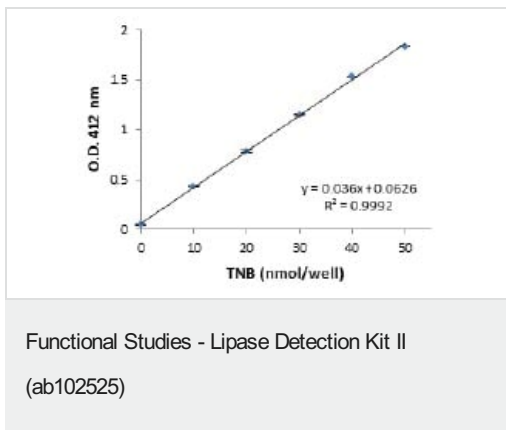
Relevance

Lipases perform essential roles in the digestion, transport and processing of dietary lipids (e.g. fats and oils) in living organisms. In humans, pancreatic lipase is the key enzyme responsible for breaking down fats in the digestive system by converting triglyceride to monoglyceride and free fatty acid. Pancreatic lipase monitoring is also used to help diagnose Crohn's disease, cystic fibrosis and celiac disease. Damage to the pancreas can exhibit a 5-10 fold increase of serum lipase levels within 24 to 48 hours.

Images



Lipase Detection in a Sample and Positive Control



TNB Standard 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.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