

Product datasheet

Lipase Assay Kit III (Fluorometric) ab118969

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

Overview

Product name	Lipase Assay Kit III (Fluorometric)
Detection method	Fluorescent
Sample type	Cell culture supernatant, Milk, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell culture media
Assay type	Enzyme activity
Sensitivity	> 0.0001 mU/well
Assay time	1h 00m
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Lipase Activity Assay Kit III (Fluorometric) (ab118969) is a lipase activity assay where lipase hydrolyzes a specific substrate to generate the methylresorufin, which can be detected fluorometrically at Ex/Em=529/600 nm. This kit can be used to detect Lipase as low as 0.1 μU/well.</p> <p>The lipase assay kit provides a rapid, simple, sensitive, and reliable test suitable for high throughput assay of Lipase activity.</p> <p>Lipase assay protocol summary:</p> <ul style="list-style-type: none">- add samples and standards to wells- add reaction mix- analyse with a microplate reader immediately and at 30-60 min
Notes	This product is manufactured by BioVision, an Abcam company and was previously called K724 Lipase Activity Fluorometric Assay Kit III. K724-100 is the same size as the 100 test size of ab118969.
Platform	Microplate reader

Properties

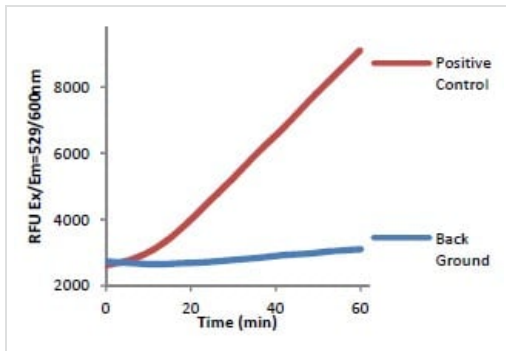
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer V	1 x 25ml
Lipase Positive Control	1 vial
Lipase Substrate I	1 x 200µl
Methylresorufin Standard	1 x 40µl

Relevance

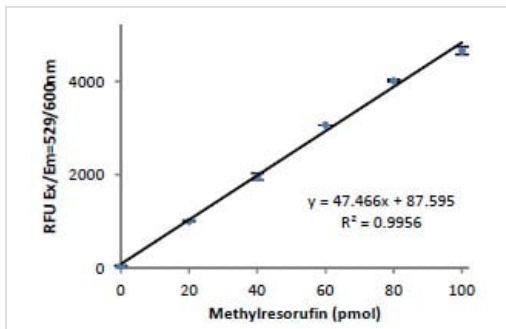
Lipases perform essential roles in the digestion, transport and processing of dietary lipids (e.g. fats and oils) in living organisms. In humans, pancreatic lipase is the key enzyme responsible for breaking down fats in the digestive system by converting triglyceride to monoglyceride and free fatty acid. 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



Positive Control and Background assayed using ab118969.

Functional Studies - Lipase Detection Kit III
(ab118969)



Example of standard curve obtained using ab118969.

Functional Studies - Lipase Detection Kit III
(ab118969)

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