

ab102525

Lipase Detection Kit II (Colorimetric)

Instructions for Use

For the rapid, sensitive and accurate measurement of Lipase activity in various samples.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

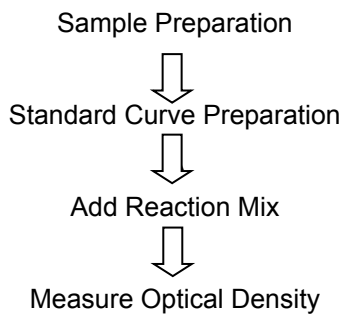
1. Overview	3
2. Protocol Summary	4
3. Components and Storage	5
4. Assay Protocol	7
5. Data Analysis	9
6. Troubleshooting	11

1. Overview

Lipases are a subclass of the esterases that catalyze the hydrolysis of ester bonds in water-insoluble, lipid substrates. Lipases perform essential roles in the digestion, transport and processing of dietary lipids (e.g. triglycerides, fats, oils) in most, if not all, living organisms. In humans, pancreatic lipases are the key enzyme responsible for breaking down fats in the digestive system by converting triglycerides to monoglycerides and free fatty acids. During the damage of the pancreas, lipase levels can rise 5 to 10-fold within 24 to 48 hours.

Abcam's Lipase Detection Kit II (Colorimetric) provides a simple, sensitive, and reliable assay for rapid analysis of Lipase in samples. In the assay, Lipases hydrolyze a specific substrate to generate a product which reacts with the DTNB to generate color ($\lambda=412$ nm). The kit is also suitable for high throughput analyses.

2. Protocol Summary



3. Components and Storage

A. Kit Components

Item	Quantity
Assay Buffer 5	25 mL
DTNB	1 vial
Lipase Substrate II	500 μ L
TNB Standard	1 vial
Lipase Positive Control	1 vial

PLEASE NOTE: Assay Buffer 5 was previously labelled as Assay Buffer V and Assay Buffer. The composition has not changed.

Store kit at -20°C , protect from light. Warm the Assay Buffer 5 to room temperature before use. Briefly centrifuge vials before opening. Read the entire protocol before performing the assay.

DTNB: Dissolve the DTNB with 1.1 ml Assay Buffer 5. Store at -20°C . Use within two months.

LIPASE SUBSTRATE II: Ready to use. Store at -20°C . Use within two months.

TNB STANDARD: Dissolve with 0.5 mL of Assay Buffer 5 to generate 5 mM TNB Standard. Stable for 2 months at -20°C.

LIPASE POSITIVE CONTROL: Dissolve the positive control with 100 µl of Assay Buffer 5. Store at -20°C. Use within two months.

B. Additional Materials Required

- Microcentrifuge
- Pipettes and pipette tips
- Colorimetric microplate reader
- 96-well plate
- Orbital shaker

4. Assay Protocol

1. Sample Preparation:

- a. **For serum samples:** Serum samples can be directly diluted in the Assay Buffer 5.
- b. **For tissue or cell samples:** Tissues or cells can be homogenized in 4 volumes of Assay Buffer 5 and centrifuged (13,000 x g, 10 min) to remove insoluble material.

Prepare test samples of up to 50 μ l/well with Assay Buffer 5 in a 96-well plate.

We suggest testing several doses of your sample to make sure readings are within the standard curve.

Note:

Mercaptans in samples will cause a high background. If the sample background is too high, the sample can be precipitated with 2 volumes of saturated ammonia sulfate. Then centrifuge, collect the precipitates and re-dissolve in the same volume of assay buffer to remove small molecule Mercaptans.

2. Positive Control (optional):

Add 5 μ l of the reconstituted Lipase Positive Control into Positive Control well and adjust the volume to 50 μ l/well with assay buffer.

3. Standard Curve Preparation:

Add 0, 2, 4, 6, 8, 10 μl of TNB Standard into a series of wells. Adjust volume to 150 μl /well with Assay Buffer 5 to generate 0, 10, 20, 30, 40, 50 nmol/well of TNB Standard.

4. Reaction Mix: Mix enough reagent for the number of assays to be performed. For each well, prepare a total 100 μl Reaction Mix.

Assay Buffer 5	85 μl
DTNB	10 μl
Lipase Substrate II	5 μl

Add 100 μl of the Reaction Mix to each well containing the Positive Controls and samples. Mix well. **(DO NOT ADD TO STANDARDS).**

5. Measurement:

Read $\text{OD}_{412\text{nm}}$ A_1 at T_1 after 3 min incubation time. Read A_2 $\text{OD}_{412\text{nm}}$ again at T_2 after incubating the reaction at 37°C for 60-90 min (or incubate longer time if the Lipase activity is low), protect from light. The OD of color generated upon formation of TNB is:

$$\Delta A_{412\text{nm}} = A_2 - A_1.$$

Note:

It is recommended to read kinetically to choose the A_1 and A_2 values which are in the linear range of the Standard Curve.

5. Data Analysis

Subtract zero Standard from all standard readings. Plot the Standard Curve.

Apply the $\Delta_{A_{412nm}}$ of samples to the standard curve to get B nmol of TNB generated in the sample reaction between T_1 and T_2 . Lipase activity in samples can then be calculated:

$$\text{Lipase Activity} = \frac{(\mathbf{B} \times \text{Dilution Factor})}{(\mathbf{T}_2 - \mathbf{T}_1) \times \mathbf{V}} = \text{nmol/min/ml} = \text{mU/ml}$$

Where:

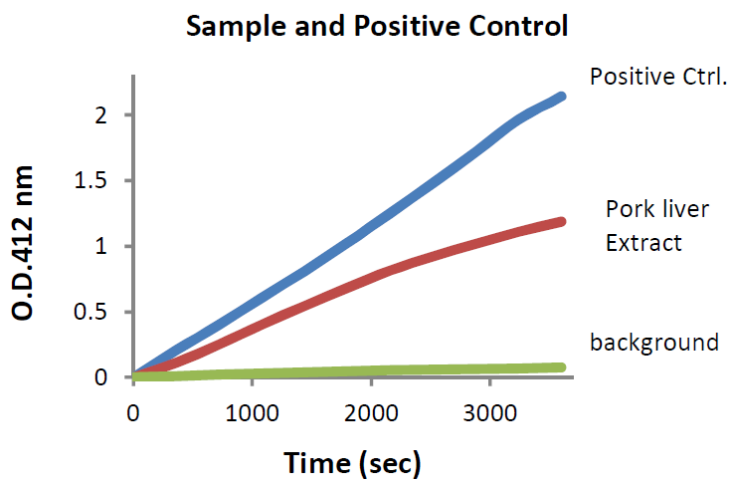
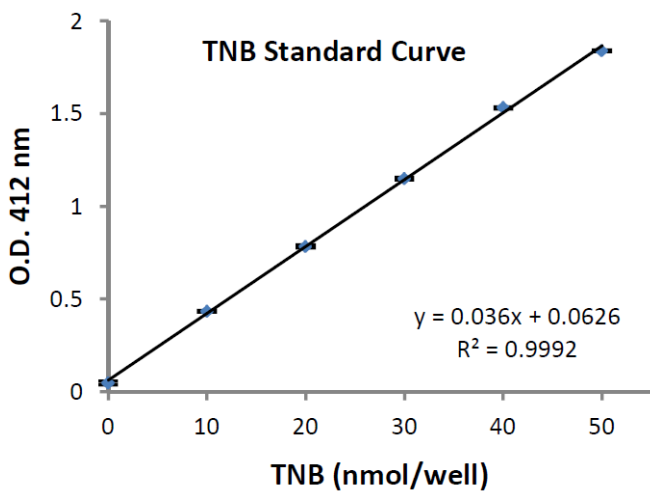
B is the TNB amount from the Standard Curve (in nmol).

T₁ is the time of the first reading (A_1) (in min).

T₂ is the time of the second reading (A_2) (in min).

V is the pre-treated sample volume (ml) added into the reaction well

Unit Definition: One unit lipase is defined as the amount of lipase which hydrolyzes the substrate and generates 1.0 μmol of TNB per minute at 37°C.



6. Troubleshooting

Problem	Reason	Solution
Assay not working	Assay buffer at wrong temperature	Assay buffer must not be chilled - needs to be at RT
	Protocol step missed	Re-read and follow the protocol exactly
	Plate read at incorrect wavelength	Ensure you are using appropriate reader and filter settings (refer to datasheet)
	Unsuitable microtiter plate for assay	Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells
Unexpected results	Measured at wrong wavelength	Use appropriate reader and filter settings described in datasheet
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Unsuitable sample type	Use recommended samples types as listed on the datasheet
	Sample readings are outside linear range	Concentrate/ dilute samples to be in linear range

Samples with inconsistent readings	Unsuitable sample type	Refer to datasheet for details about incompatible samples
	Samples prepared in the wrong buffer	Use the assay buffer provided (or refer to datasheet for instructions)
	Samples not deproteinized (if indicated on datasheet)	Use the 10kDa spin column (ab93349)
	Cell/ tissue samples not sufficiently homogenized	Increase sonication time/ number of strokes with the Dounce homogenizer
	Too many freeze-thaw cycles	Aliquot samples to reduce the number of freeze-thaw cycles
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Samples are too old or incorrectly stored	Use freshly made samples and store at recommended temperature until use
Lower/ Higher readings in samples and standards	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Out-of-date kit or incorrectly stored reagents	Always check expiry date and store kit components as recommended on the datasheet
	Reagents sitting for extended periods on ice	Try to prepare a fresh reaction mix prior to each use
	Incorrect incubation time/ temperature	Refer to datasheet for recommended incubation time and/ or temperature
	Incorrect amounts used	Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume)

Problem	Reason	Solution
Standard curve is not linear	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Pipetting errors when setting up the standard curve	Try not to pipette too small volumes
	Incorrect pipetting when preparing the reaction mix	Always prepare a master mix
	Air bubbles in wells	Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates
	Concentration of standard stock incorrect	Recheck datasheet for recommended concentrations of standard stocks
	Errors in standard curve calculations	Refer to datasheet and re-check the calculations
	Use of other reagents than those provided with the kit	Use fresh components from the same kit

For further technical questions please do not hesitate to contact us by email (technical@abcam.com) or phone (select “*contact us*” on www.abcam.com for the phone number for your region).

Technical Support

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)