

## Product datasheet

# Cholesterol/Cholesteryl Ester Assay Kit - Detection ab102515

[19 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Cholesterol/Cholesteryl Ester Assay Kit - Detection
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell Lysate, Tissue Lysate
<b>Assay type</b>	Quantitative
<b>Assay time</b>	1h 00m
<b>Product overview</b>	Cholesterol/Cholesteryl Ester Assay Kit ab102515 uses a simple method for sensitive quantification of free cholesterol, cholesteryl esters, or both using a colorimetric method.

This assay kit uses an alternative method compared to our most popular **Cholesterol/Cholesteryl Ester Assay Kit ab65359** (which relies on a cholesterol oxidase based method).

In this assay kit, free cholesterol is oxidized by cholesterol dehydrogenase to generate NADH which reacts with a sensitive probe resulting in strong absorbance at 450 nm.

The assay can detect free or total cholesterol depending upon whether cholesterol esterase is used to hydrolyze cholesterol esters present.

Cholesteryl ester can be determined by subtracting the value of free cholesterol from the total cholesterol (cholesterol plus cholesteryl esters).

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called K623 Total Cholesterol and Cholesteryl Ester Colorimetric Assay Kit II. K623-100 is the same size as the 100 test size of ab102515.

Cholesterol is an essential molecule in all animal life. It has been involved in both normal development and diseases.

**Platform** Microplate reader

### Properties

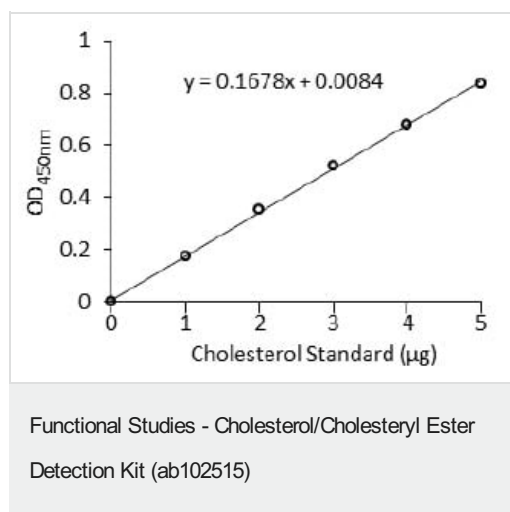
**Storage instructions** Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer II	1 x 25ml
Cholesterol Enzyme Mix	1 vial
Cholesterol Esterase	1 vial
Cholesterol Standard	1 x 100µl
Developer Solution III	1 vial

## Relevance

Cholesterol is an essential molecule in all animal life. It is involved in both normal development and diseases. A high level of serum cholesterol is an indicator for diseases such as heart disease. Cholesteryl ester is an ester of cholesterol. The ester bond is formed between the carboxylate group of a fatty acid and the hydroxyl group of cholesterol.

## Images



Cholesterol Standard Curve: Cholesterol/Cholesteryl Ester was quantified according to the kit instructions. Background from the 0 standard reading (without cholesterol) has been subtracted from all readings.

**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