

Product datasheet

Mouse free cholesterol ELISA Kit ab285294

[1 Image](#)

Overview

Product name	Mouse free cholesterol ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma, Tissue Homogenate
Assay type	Sandwich (quantitative)
Range	1.25 ng/ml - 80 ng/ml
Assay time	0h 90m
Assay duration	Multiple steps standard assay
Product overview	Mouse free cholesterol ELISA kit ab285294 (previously called QuickDetect free cholesterol (fCHOL) ELISA Kit, K4438) is a sandwich ELISA assay for measuring the quantity of free Cholesterol in mouse serum, plasma, cell culture supernatants cell culture, tissue homogenates and other biological fluids. The entire process can be performed within 90 minutes with only a single incubation and a wash step respectively, resulting in fewer handling steps, fewer errors and more consistent results.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called K4438 QuickDetect™ free cholesterol (fCHOL) (Mouse) ELISA Kit. K4438-100 is the same size as the 96 test size of ab285294.
Platform	Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions Store at +4°C. Please refer to protocols.

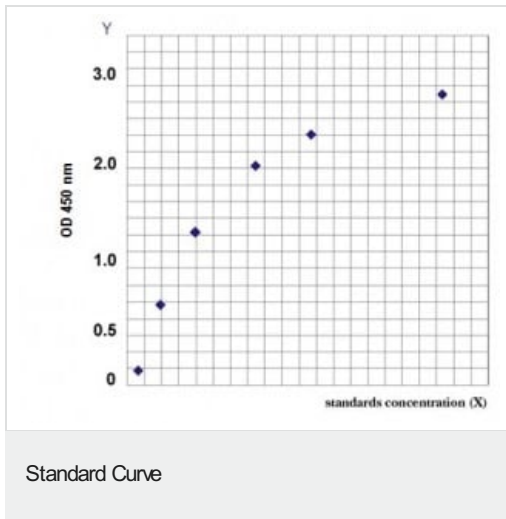
Components	96 tests	96 tests
20X Wash buffer	1 x 20ml	1 x 20ml
Chomogen Solution A	1 x 6ml	1 x 6ml
Chomogen Solution B	1 x 6ml	1 x 6ml
HRP-Conjugate Reagent	1 x 6ml	1 x 6ml

Components	96 tests	96 tests
Micro ELISA Plate	1 unit	1 unit
Microplate Sealer	2 units	2 units
Special diluent	1 x 6ml	1 x 6ml
Standard (80 ng/ml)	1 x 0.6ml	1 x 0.6ml
Standard diluent	1 x 6ml	1 x 6ml
Stop Solution	1 x 6ml	1 x 6ml

Relevance

Cholesterol is a waxy substance that occurs naturally in all parts of the body and that the body needs to function normally. It is present in cell walls or membranes everywhere in the body, including the brain, nerves, muscle, skin, liver, intestines, and heart. The body uses cholesterol to produce many hormones, vitamin D, and the bile acids that help to digest fat. It takes only a small amount of cholesterol in the blood to meet these needs. However, if you have too much cholesterol in your bloodstream, it can lead to atherosclerosis, a condition in which fat and cholesterol are deposited in the walls of the arteries in many parts of the body, including the coronary arteries feeding the heart. In time, narrowing of the coronary arteries by atherosclerosis can produce the signs and symptoms of heart disease, including angina and heart attack.

Images



Typical Standard Curve: These standard curves are for demonstration only. A standard curve must be run with each assay.

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