

Triglyceride Assay Kit (Fluorometric, Reducing Samples) ab178780

★★★★★ [1 Abreviews](#) [11 References](#) [2 图像](#)

概述

产品名称	Triglyceride Assay试剂盒(Fluorometric, Reducing Samples)
检测方法	Fluorescent
样品类型	Saliva, Serum, Plasma, Tissue, Adherent cells, Suspension cells
检测类型	Quantitative
灵敏度	< 0.4 μ M
检测时间	1h 00m
种属反应性	与反应: Mammals, Other species
产品概述	<p>Triglyceride Assay Kit (ab178780) is a sensitive, easy assay to measure triglyceride concentration in mammalian samples. In this triglyceride assay protocol, triglycerides are hydrolyzed to glycerol and fatty acid. The glycerol reacts with Triglyceride Enzyme Mix to form an intermediate product which in turn reacts with the probe and developer to generate fluorescence (Ed/Em 535/587 nm). The generated fluorescence is directly proportional to the amount of triglycerides.</p>

Triglyceride Assay Kit (Fluorometric, Reducing Samples) ab178780 is suitable for measuring triglyceride levels in samples which contain reducing substances that may interfere with oxidase-based assays such as [Triglyceride Assay Kit \(ab65336\)](#).

Triglyceride assay protocol summary:

- add samples and standards to wells
- add lipase and incubate for 20 min
- add reaction mix and incubate for 30 min
- analyze with a microplate reader.

Please note: the general range is 0-10 nmol (colorimetric) and 0-1 nmol (fluorometric).

Target background

Triglycerides are the main constituent of vegetable oil, animal fat, LDL and VLDL, and play an important role as transporters of fatty acids as well as serving as an energy source. Triglycerides are broken down into fatty acids and glycerol, after which both can serve as substrates for energy

producing and metabolic pathways. High blood levels of triglycerides are implicated in atherosclerosis, heart disease and stroke as well as in pancreatitis.

Review our [Metabolism Assay Guide](#) to learn about assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also about how to assay metabolic function in live cells using your plate reader.

说明

This product is manufactured by BioVision, an Abcam company and was previously called K614 PicoProbe™ Triglyceride (TG) Fluorometric Assay Kit. K614-100 is the same size as the 100 test size of ab178780.

How other researchers have used Triglyceride Assay Kit ab178780

The Triglyceride assay kit has been used in publications in a variety of sample types, including:

- Human: Hepatocyte cell/cell line extract^{1,2}
- Mouse: Plasma³, Skeletal muscle extract⁴, Serum⁵

References: 1 - [Liu Q et al. 2021](#), 2 - [Hu LK et al. 2021](#), 3 - [Panajatovic MV et al.](#), 4 - [Panajatovic MV et al.](#), 5 - [Maier JI et al. 2021](#)

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

平台

Microplate reader

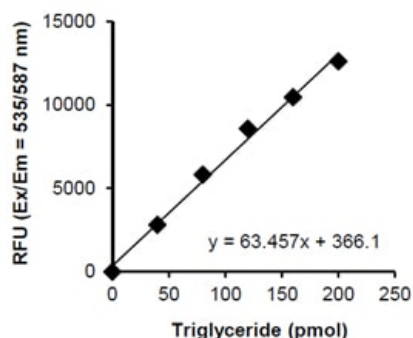
性能

存放说明

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

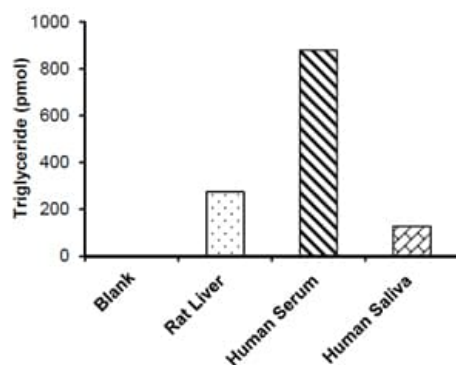
组件	100 tests
Assay Buffer V	1 x 25ml
Cholesterol Esterase	1 vial
PicoProbe I	1 x 0.4ml
Substrate Mix I	1 vial
Triglyceride Enzyme Mix	1 vial
Triglyceride Standard	1 x 0.3ml

图片



Example of triglyceride standard curve following assay protocol. Please note that this is an example curve and should not be used for experimental purposes.

Triglyceride Standard Curve



Mesurament of TG levels in rat liver (~15 µg), human serum (1 µL) and saliva (10 µL). Assays were performed according to kit protocol.

Measurement of triglyceride levels in samples

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors