

Product datasheet

Free Glycerol Assay Kit (Fluorometric, High Sensitivity)
ab174092

4 Images

Overview

Product name	Free Glycerol Assay Kit (Fluorometric, High Sensitivity)
Detection method	Fluorescent
Sample type	Serum, Plasma, Other biological fluids, Tissue, Adherent cells, Suspension cells
Assay type	Quantitative
Sensitivity	0.04 nmol/well
Species reactivity	Reacts with: Other species, Mammals
Product overview	Abcam's PicoProbe Free Glycerol Assay Kit (ab174092) is simple, sensitive and easy to use tool to detect small amounts of glycerol.

This assay is suitable for measuring trace amount of glycerol in samples containing reducing substances, which may interfere with oxidase-based assays. In the assay, glycerol reacts with the enzyme mix to form an intermediate, which is subsequently oxidized with the production of fluorescence. The fluorescence intensity is directly proportional to the amount of glycerol.

This assay kit can detect glycerol amount less than 40 pmol.

Visit our [FAQs](#) page for tips and troubleshooting.

Notes Glycerol is a central component for synthesis of all lipids; it acts as a backbone for triglycerides and phospholipids, which plays an important role for cell membrane's structure. Due to its low toxicity, glycerol is widely used in pharmaceutical, food and cosmetic industries.

Platform Microplate reader

Properties

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

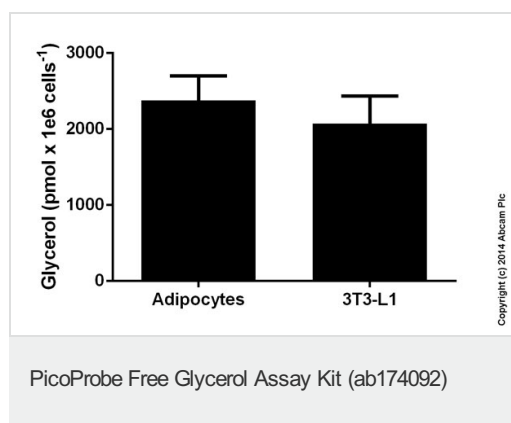
Components	Identifier	100 tests
Glycerol Assay Buffer	WM	1 x 25ml
Glycerol Developer		1 vial

Components	Identifier	100 tests
Glycerol Enzyme Mix (lyophilized)	Green	1 vial
Glycerol Standard	Yellow	1 x 0.2ml
PicoProbe	Blue	1 x 0.4ml

Relevance

Glycerol is widely used in foods, beverages, solvents, pharmaceutical and cosmetic products, etc. There is broad interest in quantification of glycerol for research and development.

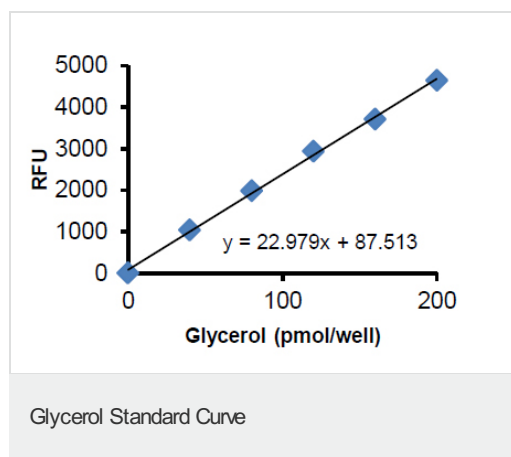
Images



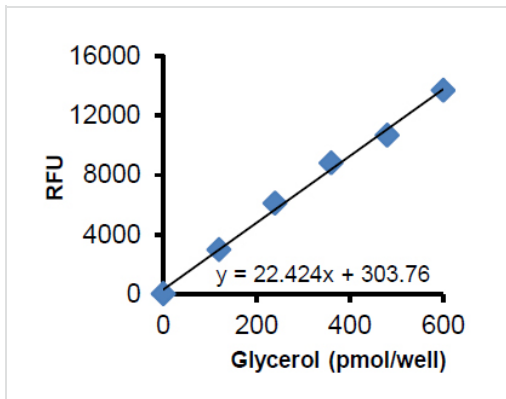
Free Glycerol measured in cell lysates showing quantity (pmol) per 1 mln cells.

Samples with the concentration of 1 e7 cells/mL were used.

Samples were diluted 2-6 fold.

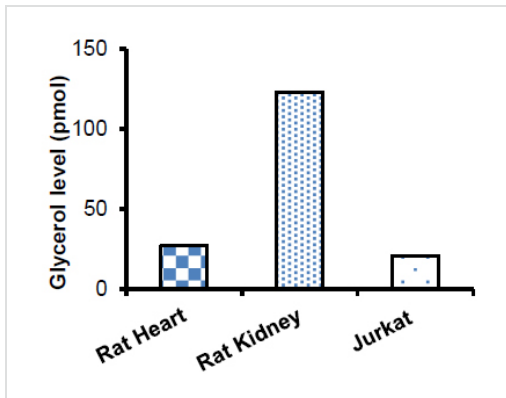


Example of Glycerol standard curve following assay's protocol.



Example of Glycerol standard curve following assay's protocol.

Glycerol Standard Curve



Measurement of glycerol levels in rat heart (5 µg protein), rat kidney (5 µg) and Jurkat cell lysate (10 µg) sample following PicoProbe Free Glycerol Assay Kit (ab174092) protocol. This is example data only.

Measurement of Glycerol Levels

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