

Product datasheet

Free Glycerol Assay Kit ab65337

[29 References](#) [5 Images](#)

Overview

Product name	Free Glycerol Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Cell culture supernatant, Serum, Plasma, Other biological fluids, Tissue Extracts
Assay type	Quantitative
Sensitivity	> 1 μ M
Range	1 μ M - 10000 μ M
Assay time	0h 40m
Product overview	Glycerol Assay Kit (ab65337) provides a sensitive, easy assay to measure free glycerol concentration.

In the glycerol assay protocol, glycerol is enzymatically oxidized to generate a product which reacts with the probe to generate color ($\lambda = 570$ nm) and fluorescence (Ex/Em = 535/587 nm).

This glycerol assay kit can detect 50 pmol-10 nmol (or ~1-10000 μ M range) of glycerol.

Glycerol assay protocol summary:

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

Notes This product is manufactured by BioVision, an Abcam company and was previously called K630 Free Glycerol Colorimetric/Fluorometric Assay Kit. K630-100 is the same size as the 100 test size of ab65337.

Platform Microplate reader

Properties

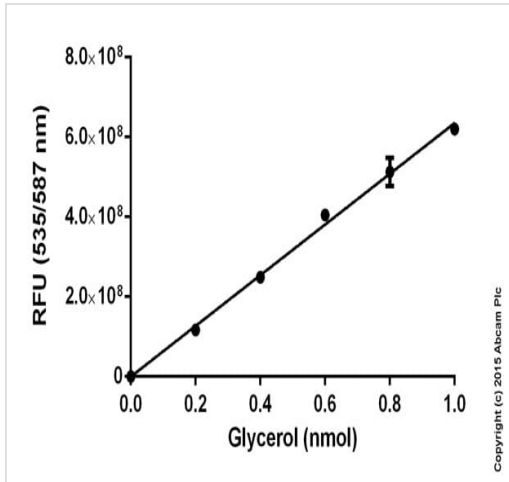
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer V	1 x 25ml

Components	100 tests
Enzyme Mix VI	1 vial
Glycerol Standard	1 x 200µl
OxiRed Probe	1 x 200µl

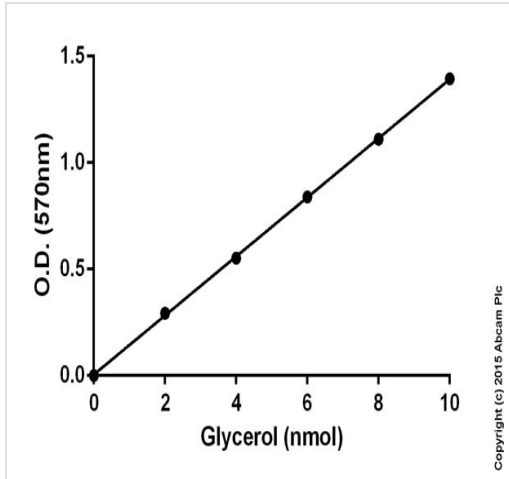
Relevance Glycerol is widely used in foods, beverages, solvents, pharmaceutical and cosmetic products, etc. There is broad interest in quantification of glycerol for research and development.

Images



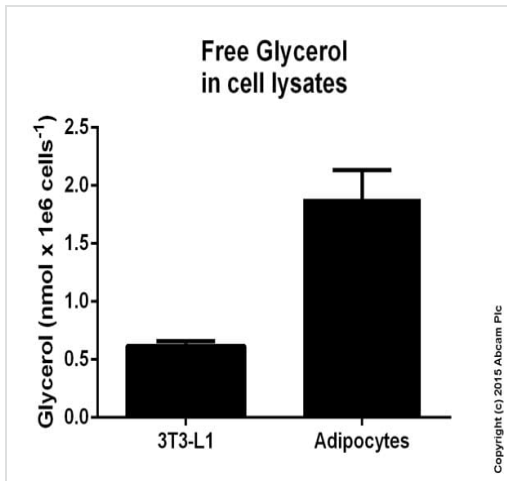
Standard curve: mean of duplicates (+/- SD) with background reads subtracted

Functional Studies - Free Glycerol Assay Kit (ab65337)



Standard curve: mean of duplicates (+/- SD) with background reads subtracted

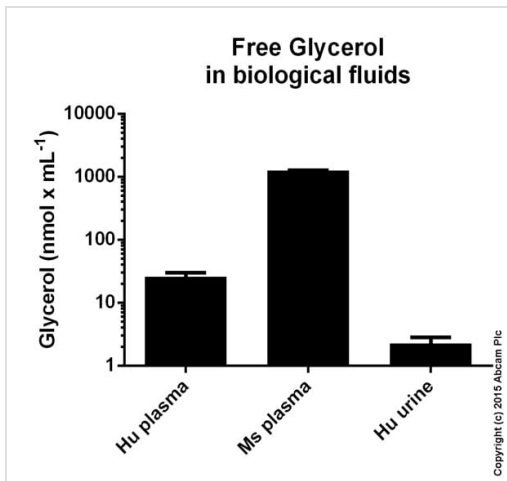
Functional Studies - Free Glycerol Assay Kit (ab65337)



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

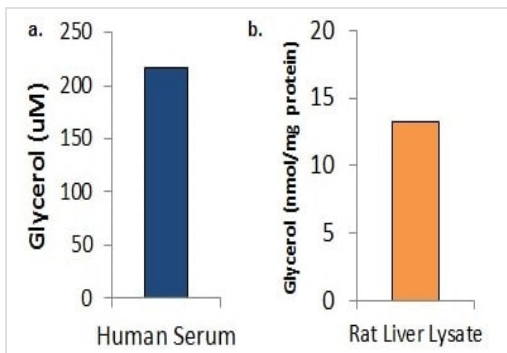
Samples with the concentration of 1e7 cells/mL were used. Samples were diluted 2-8 fold and measured fluorometrically.

Functional Studies - Free Glycerol Assay Kit (ab65337)



Free Glycerol measured in biological fluids showing quantity (nmol) per mL of tested sample. Samples were diluted 2-8 fold and measured colorimetrically.

Functional Studies - Free Glycerol Assay Kit (ab65337)



a. Measurement of free glycerol in pooled human serum (10 µL). b. Measurement of free glycerol in rat liver lysate (500 µg).

Free Glycerol Assay Kit - Functional Studies

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