

### Free Glycerol Assay Kit ab65337

**29 References**    **画像数 5**

#### 製品の概要

製品名	Free Glycerol Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell culture supernatant, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
検出感度	> 1 µM
検出範囲	1 µM - 10000 µM
全工程の試験時間	0h 40m
製品の概要	<p>Glycerol Assay Kit (ab65337) provides a sensitive, easy assay to measure free glycerol concentration.</p> <p>In the glycerol assay protocol, glycerol is enzymatically oxidized to generate a product which reacts with the probe to generate color (<math>\lambda = 570</math> nm) and fluorescence (Ex/Em = 535/587 nm).</p> <p>This glycerol assay kit can detect 50 pmol-10 nmol (or ~1-10000 µM range) of glycerol.</p> <p>Glycerol assay protocol summary:</p> <ul style="list-style-type: none"> <li>- add samples and standards to wells</li> <li>- add reaction mix and incubate for 30 min</li> <li>- analyze with a microplate reader</li> </ul>
特記事項	<p>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.</p> <p>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.</p> <p>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.</p>
試験プラットフォーム	Microplate reader

#### 製品の特性

保存方法

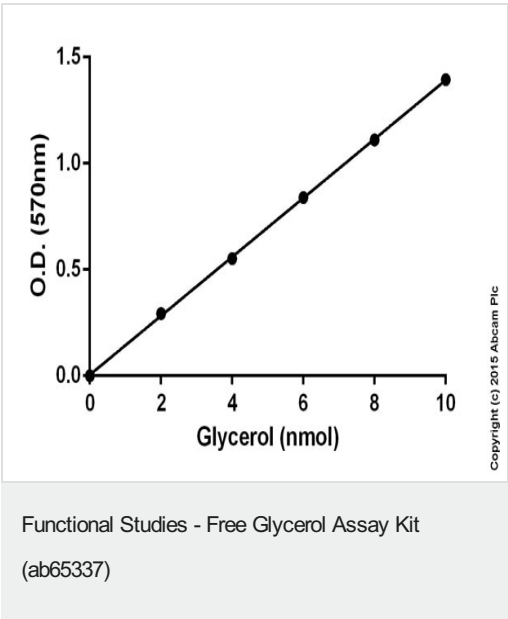
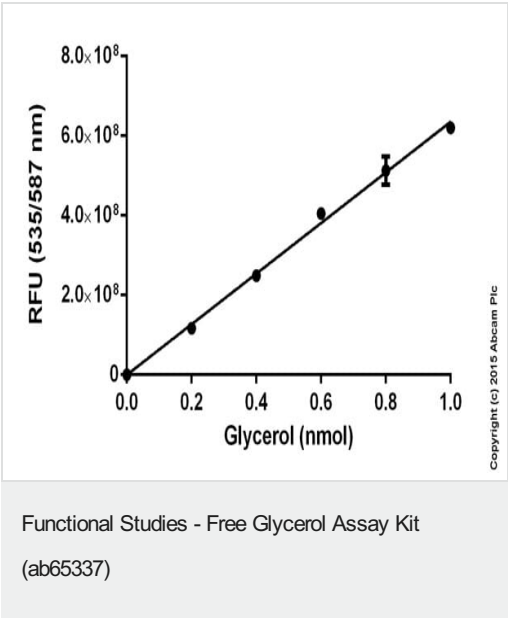
Store at -20°C. Please refer to protocols.

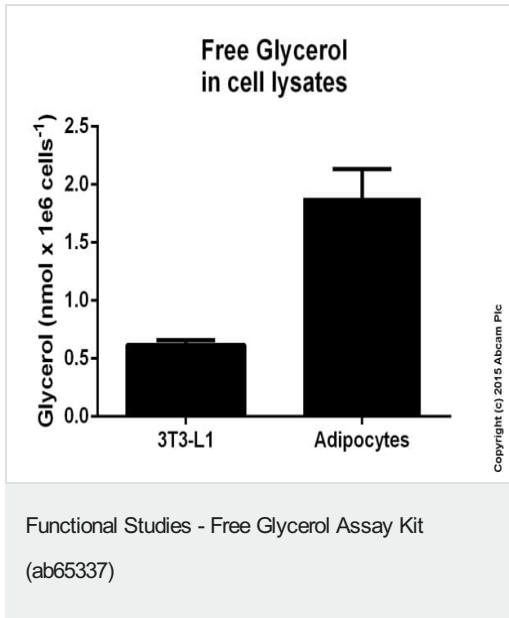
内容	100 tests
Assay Buffer V	1 x 25ml
Enzyme Mix VI	1 vial
Glycerol Standard	1 x 200µl
OxiRed Probe	1 x 200µl

関連性

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.

画像

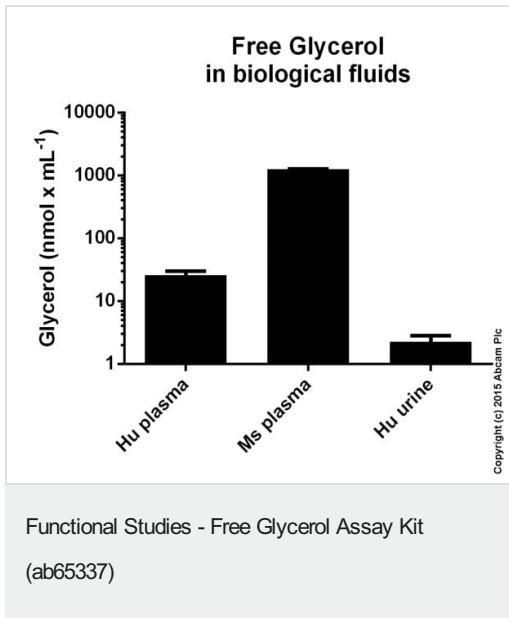




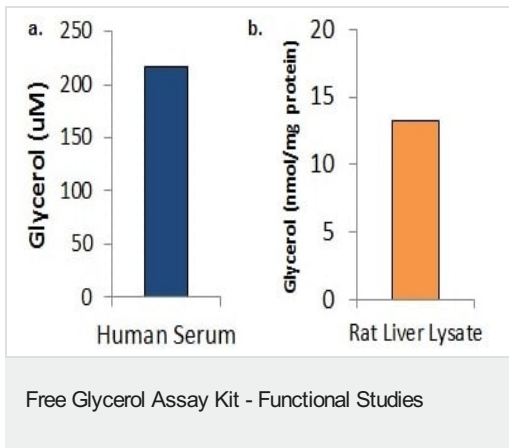
Free Glycerol measured in cell culture lysates showing quantity (nmol) per 1 mln cells.

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

Samples were diluted 2-8 fold and measured fluorometrically.



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



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

**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.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

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