

Free Glycerol Assay Kit II (Colorimetric) ab155899

[1 References](#) [画像数 1](#)

製品の概要

製品名	Free Glycerol Assay Kit II (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Tissue, Adherent cells, Suspension cells, Tissue Homogenate
アッセイタイプ	Quantitative
検出感度	< 20 μ M
検出範囲	2 nmol/well - 10 nmol/well
全工程の試験時間	1h 30m
種交差性	交差種: Mammals, Other species
製品の概要	Abcam's Free Glycerol Assay kit II (ab155899) is suitable for measuring free glycerol levels in samples that contain reducing substances, which may interfere with oxidase-based assays. In this assay, Glycerol in the presence of Glycerol Enzyme Mix is converted to an intermediate, which reduces a colorless Probe to a colored product with strong absorbance at 450 nm. Free Glycerol Assay Kit II is simple, rapid and high-throughput adaptable. This assay kit can detect less than 20 μ M of free Glycerol in various biological samples.

Detection method: colorimetric – based on enzyme activity

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K634 Free Glycerol Colorimetric Assay Kit II. K634-100 is the same size as the 100 test size of ab155899.</p> <p>Glycerol is a central component of most lipids. It acts as a backbone for triglycerides and phospholipids, which play an important role in metabolism and cell membrane structure. Due to its low toxicity, glycerol is widely used in the pharmaceutical, food and cosmetic industries.</p>
------	--

試験プラットフォーム	Microplate reader
------------	-------------------

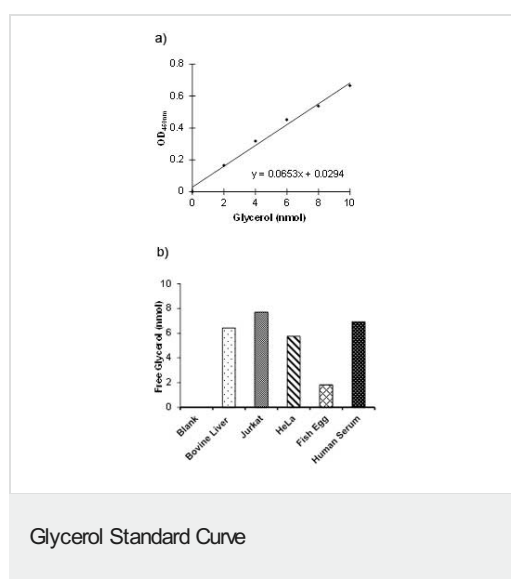
製品の特性

保存方法	Store at -20°C. Please refer to protocols.
------	--

内容	100 tests
Developer Solution III	1 vial
Glycerol Assay Buffer	1 x 25ml
Glycerol Enzymatic Developer	1 vial
Glycerol Standard	1 x 0.2ml

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



Glycerol Standard Curve [a]. Measurement of free glycerol in bovine liver (50 μ g), Jurkat cells (40 μ g), HeLa cells (30 μ g), fish egg (58 μ g) & human serum (2 μ l) [b]. Assays were performed following kit protocol.

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