

Lipolysis Assay Kit (Fluorometric) ab185434

1 References 画像数 3

製品の概要

製品名	Lipolysis Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Cell culture supernatant
アッセイタイプ	Quantitative
検出感度	< 0.02 nmol/well
種交差性	交差種: Mammals, Other species
製品の概要	<p>Lipolysis Assay Kit (Fluorometric) ab185434 is simple, easy-to-use and highly sensitive.</p> <p>This lipolysis assay measures glycerol released from 3T3-L1 cells as early as 1 hour after induction of lipolysis. It is suitable for measuring trace amounts of glycerol from samples. In the assay, glycerol reacts with Glycerol Enzyme Mix to form an intermediate that in turn reacts with Glycerol Developer and PicoProbe™ to generate the fluorescence product. The fluorescence intensity is directly proportional to the amount of glycerol.</p> <p>The lipolysis assay kit can detect less than 20 pmol of glycerol.</p>
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K578 PicoProbe™ Lipolysis (3T3-L1) Fluorometric Assay Kit. K578-100 is the same size as the 100 test size of ab185434.</p> <p>Lipolysis is the hydrolysis of triglycerides within the cell into glycerol and free fatty acids. The glycerol and free fatty acids are then released into the bloodstream or culture media. Lipolysis occurs in essentially all cells, but is most abundant in white and brown adipose tissue.</p>
試験プラットフォーム	Microplate reader

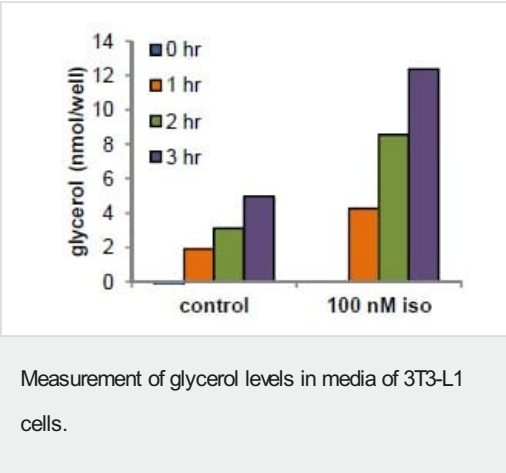
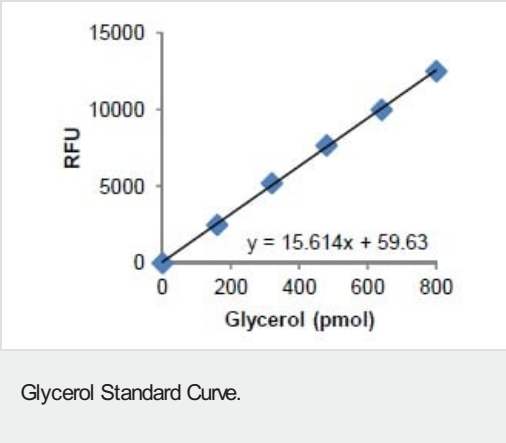
製品の特性

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

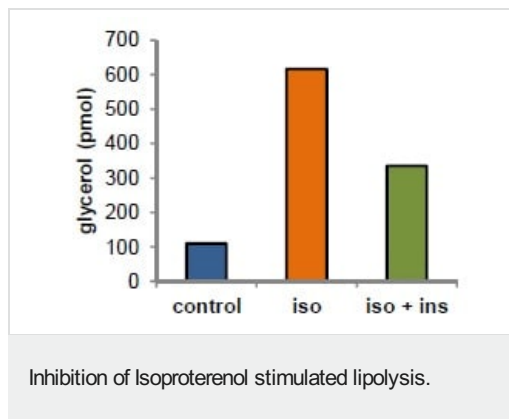
内容	100 tests
Glycerol Assay Buffer	1 x 25ml

内容	100 tests
Glycerol Developer	1 vial
Glycerol Enzyme Mix	1 vial
Glycerol Standard (100 mM)	1 x 0.2ml
Isoproterenol (10mM)	1 x 50μl
Lipolysis Assay Buffer	1 x 17ml
Lipolysis Wash Buffer	1 x 22ml
PicoProbe	1 x 0.4ml

画像



5 μl of 3T3-L1 cells treated with vehicle control (H₂O) or 100 nM Isoproterenol for 0-3 hours. Measurements for 0 hours were undetectable.



Inhibition of Isoproterenol (100 nM) stimulated lipolysis by treatment with 100 nM insulin (measured using 5 µl of media of 3T3-L1 cells). Control was vehicle (H₂O) treated cells.

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