

Lipolysis Assay Kit (Colorimetric) ab185433

★★★★★ [1 Abreviews](#) [14 References](#) [画像数 2](#)

製品の概要

製品名	Lipolysis Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant
アッセイタイプ	Quantitative
検出感度	< 0.2 nmol/well
検出範囲	2 nmol/well - 10 nmol/well
全工程の試験時間	0h 40m
種交差性	交差種: Mammals, Other species
製品の概要	<p>Abcam's Lipolysis Assay Kit (Colorimetric) (ab185433) is simple and easy-to-use. The assay measures glycerol released from 3T3-L1 cells after induction of lipolysis using colorimetric method. The color intensity is directly proportional to the amount of glycerol. This assay kit can detect less than 200 pmol of glycerol.</p>

Lipolysis assay protocol summary:

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

特記事項 This product is manufactured by BioVision, an Abcam company and was previously called K577 Lipolysis (3T3-L1) Colorimetric Assay Kit. K577-100 is the same size as the 100 test size of ab185433.

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. Deficiencies in lipolysis lead to increased intracellular lipid accumulation, resulting in abnormal cellular physiology, hyperlipidemia, and insulin resistance. Lipolysis can be induced by catecholamine and certain hormones. The kit includes synthetic catecholamine, Isoproterenol, which activates β -adrenergic receptors. This leads to activation of adenylate cyclase, which catalyzes the conversion of ATP to cAMP. cAMP then serves as a second messenger to activate hormone-sensitive lipase, which hydrolyzes the triglycerides. This pathway can be inhibited by insulin.

試験プラットフォーム Microplate reader

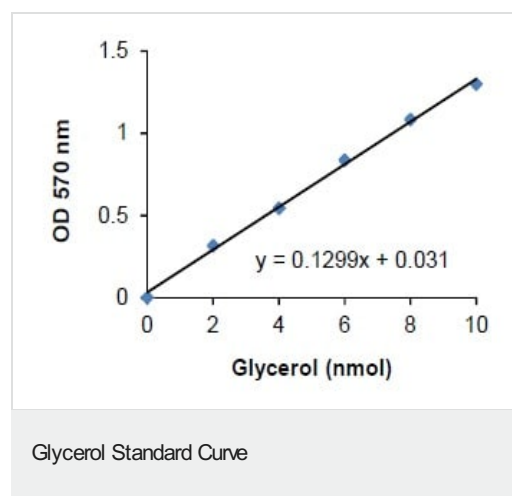
製品の特性

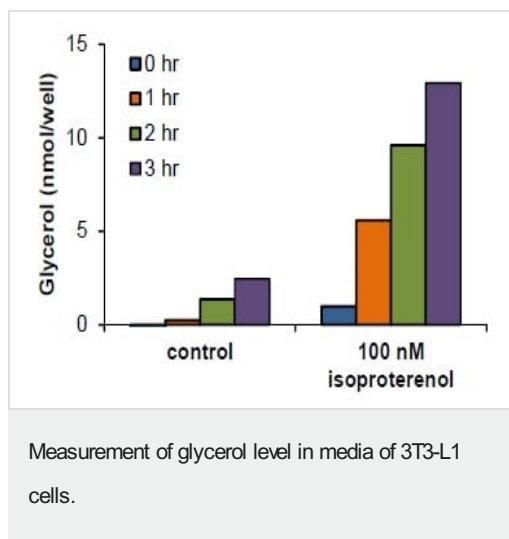
保存方法

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 0.2ml
Isoproterenol	1 x 50µl
Buffer IV	1 x 17ml
Wash Buffer VIII	1 x 22ml
OxiRed Probe	1 x 0.2ml

画像





50 μ l of 3T3-L1 cells treated with vehicle control (H_2O) or 100 nM Isoproterenol for 0-3 hours.

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