

LDH-Cytotoxicity Assay Kit (Fluorometric) ab197004

[10 References](#) [3 图像](#)

概述

| | |
|-------|--|
| 产品名称 | LDH-Cytotoxicity Assay试剂盒(Fluorometric) |
| 检测方法 | Fluorescent |
| 样品类型 | Adherent cells, Suspension cells |
| 检测类型 | Enzyme activity (quantitative) |
| 灵敏度 | 100 cells/well |
| 检测时间 | 0h 20m |
| 种属反应性 | 与反应: Mammals, Other species |
| 产品概述 | LDH Assay Kit (Cytotoxicity, Fluorometric) (ab197004) provides a sensitive, quick, and easy way to detect LDH (lactate dehydrogenase) released from damaged cells. |

In the LDH assay, LDH converts lactate to pyruvate and NADH, which reduces a proprietary probe to an intensely fluorescent product (Ex/Em = 535/587 nm).

The amount of fluorescence is directly proportional to the number of damaged cells.

The assay is adaptable to high-throughput format and can be completed in less than 20 min. Sensitivity: ~ 100 cells. 5µl of cell culture media required per assay.

说明 This product is manufactured by BioVision, an Abcam company and was previously called K314 PicoProbe™ LDH-Cytotoxicity Fluorometric Assay Kit. K314-500 is the same size as the 500 test size of ab197004.

Alternative LDH assays

If you would like to use a fluorometric assay, please refer to [LDH-Cytotoxicity Assay Kit \(Fluorometric\) \(ab197004\)](#).

This kit is more sensitive than the [colorimetric LDH Cytotoxicity Assay Kit ab65391](#).

To measure LDH activity in sample types such as serum, plasma, and cell lysates, we recommend [LDH assay kit ab102526](#).

Related products and guides to cytotoxicity / cell viability / proliferation assays

Review our [cell health assay guide](#) to learn about our other kits to perform a [cell viability assay](#), [cytotoxicity assay](#) or [cell proliferation assay](#).

平台 Microplate reader

性能

存放说明Store at -20°C. Please refer to protocols.

| 组件 | 500 tests | 500 tests |
|----------------------|-----------|-----------|
| Developer Mix I | 1 vial | 1 vial |
| LDH Assay Buffer | 1 x 50ml | 1 x 50ml |
| LDH Positive Control | 1 x 20µl | 1 x 20µl |
| Lysis Buffer II | 1 x 5ml | 1 x 5ml |
| PicoProbe III | 1 x 2ml | 1 x 2ml |

通路Fermentation; pyruvate fermentation to lactate; (S)-lactate from pyruvate: step 1/1.

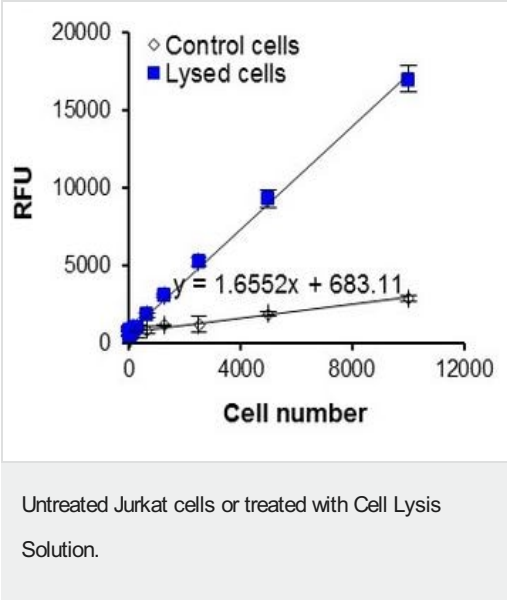
疾病相关Defects in LDHA are the cause of glycogen storage disease type 11 (GSD11) [MIM:612933]. A metabolic disorder that results in exertional myoglobinuria, pain, cramps and easy fatigue.

序列相似性Belongs to the LDH/MDH superfamily. LDH family.

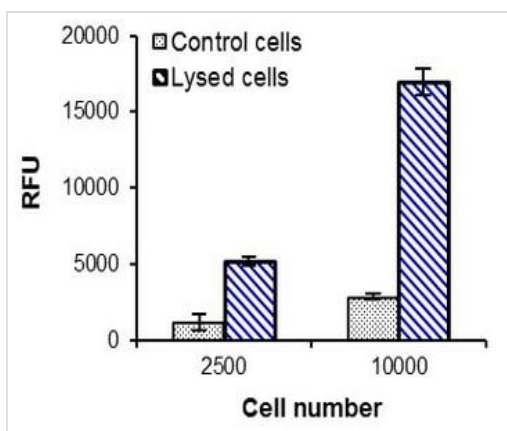
翻译后修饰ISGylated.

细胞定位Cytoplasm.

图片

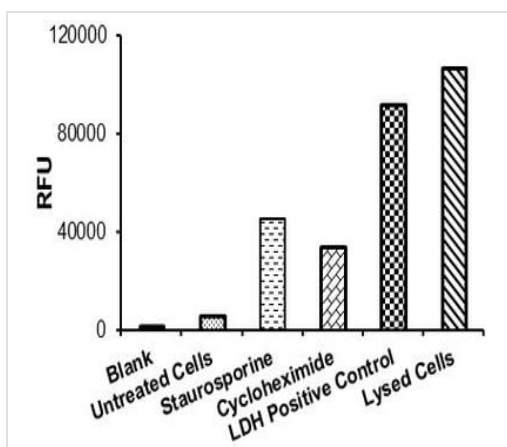


Untreated Jurkat cells (low control) or treated with Cell Lysis Solution for 30 min (high control).



Relative fluorescence units of untreated Jurkat cells (low control) or treated with Cell Lysis Solution for 30 min (high control).

Relative Fluorescence Units of untreated Jurkat cells or treated with Cell Lysis Solution.



Overnight treatment of HeLa cells with 100 μ M of staurosporine or 3 μ M of cycloheximide. LDH released into the medium was measured along with blank, untreated cells (low control), LDH Positive Control, and lysed cells (high control)

Overnight treatment of HeLa cells with staurosporine or cycloheximide.

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.cn/abpromise> or contact our technical team.

Terms and conditions

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