

LDH Assay Kit (Cytotoxicity) ab65391

35 References **2 图像**

概述

产品名称	LDH Assay试剂盒 (Cytotoxicity)
样品类型	Adherent cells, Suspension cells
检测类型	Enzyme activity
检测时间	1h 00m
产品概述	For a more sensitive LDH Assay Kit to use as a cytotoxicity assay, we recommend ab65393 .

LDH Assay Kit (Cytotoxicity) ab65391 provides a fast and simple method for quantitating cytotoxicity based on the measurement of activity of lactate dehydrogenase (LDH) released from damaged cells.

The assay takes ~0.5-1 hr. LDH activity can be determined by a coupled enzymatic reaction: LDH oxidizes lactate to pyruvate which then reacts with tetrazolium salt INT to form formazan. The increase in the amount of formazan produced in culture supernatant directly correlates to the increase in the number of lysed cells. The formazan dye is water-soluble and can be detected by spectrophotometer at 500 nm.

This kit was previously called LDH-Cytotoxicity Assay Kit I.

说明 This product is manufactured by BioVision, an Abcam company and was previously called K311 LDH-Cytotoxicity Colorimetric Assay Kit. K311-400 is the same size as the 400 test size of ab65391.

Cell death or cytotoxicity is classically evaluated by the quantification of plasma membrane damage. Unlike many other cytoplasmic enzymes which exist in many cells either in low amount (e.g., alkaline and acid phosphatase) or unstable, LDH is a stable cytoplasmic enzyme present in all cells and rapidly released into the cell culture supernatant upon damage of the plasma membrane.

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

平台 Microplate reader

性能

存放说明

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

组件	400 tests	8000 tests
LDH Catalyst Reagent	1 vial	20 vials
Dye Solution I	1 x 45ml	20 x 45ml

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

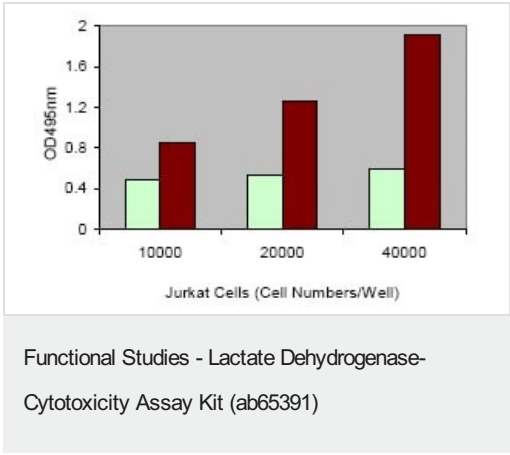
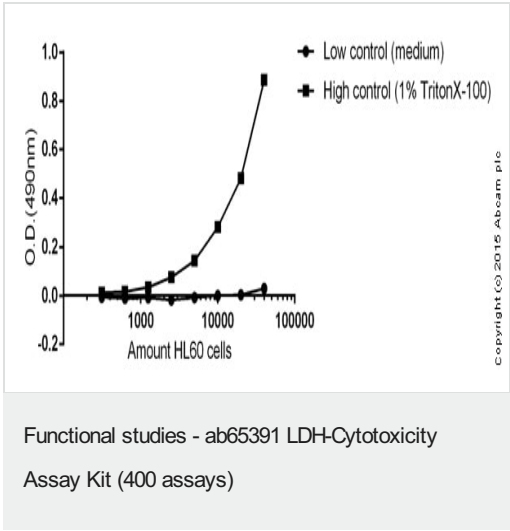
疾病相关 Defects in LDHA are the cause of glycogen storage disease type 11 (GSD11) [MIM:612933]. A metabolic disorder that results in exertional myoglobinuria, pain, cramps and easy fatigue.

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



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