

# ALDH Activity Assay Kit (Colorimetric) ab155893

★★★★★ [1 Abreviews](#) [11 References](#) [画像数 3](#)

## 製品の概要

製品名	ALDH Activity Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma, Cell Lysate, Tissue Culture Media
アッセイタイプ	Enzyme activity (quantitative)
検出感度	< 0.5 mU/well
全工程の試験時間	0h 30m
種交差性	交差種: Mammals, Other species
製品の概要	ALDH Activity Assay Kit (Colorimetric) (ab155893) is a simple, fast and reliable method to quantify the ALDH enzymatic activity. In this assay, acetaldehyde is oxidized by ALDH generating NADH which then reduces a colorless probe to a colored product with strong absorbance at 450 nm. The assay can detect < 0.5 mU of ALDH activity in a variety of samples.

ALDH assay protocol summary:

- add samples and standards to wells
- add reaction mix and incubate for 20-60 min
- analyze with a microplate reader at 5 min and 20-60 min

## 特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K731 Aldehyde Dehydrogenase Activity Colorimetric Assay Kit. K731-100 is the same size as the 100 test size of ab155893.

The NAD-dependent Aldehyde Dehydrogenase (ALDH) plays a vital role in cellular detoxification. It oxidizes various aldehydes and generates the corresponding carboxylic acid. ALDH have been found in every cellular compartment. Based on its structure and function, ALDH comprises 3 major classes in mammals: Class 1 and Class 3 (the tumor form) are located in the cytosol and include both constitutive and induced forms; Class 2 is located in the mitochondria and only exists as the constitutive form. In humans, the ALDH superfamily consists of 19 genes. The mutation of ALDH genes (loss of function) causes human diseases such as Type II hyperprolinemia, pyridoxine-dependent seizure and hyperammonemia. Recent studies show that increased ALDH activity leads to several types of malignancies, serves as a cancer stem cell marker and correlates with poor prognosis.

## 試験プラットフォーム

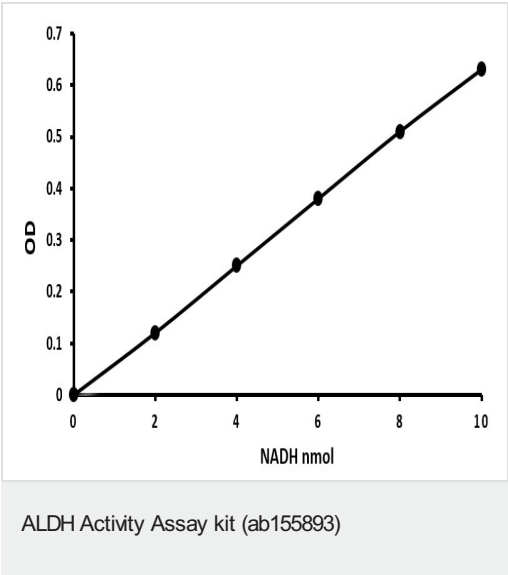
Microplate reader

## 製品の特性

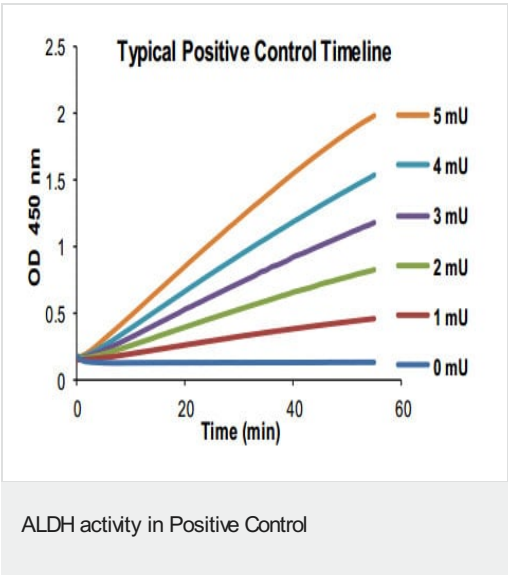
保存方法Store at -20°C. Please refer to protocols.

内容	100 tests
Acetaldehyde	1 x 0.5ml
Assay Buffer XXVIII	1 x 25ml
ALDH Positive Control	1 vial
Developer Solution III	1 vial
NADH Standard I	1 vial

画像



Standard curve (60 minutes incubation) showing mean of duplicates (+/- SD) with background reads subtracted.



Typical OD of the diffrent positive contols mesuared over 1 hour.



ALDH activity measured in human biological fluids ( $T_1=5$  min;  $T_2=60$  min), showing activity (mU) per mL of tested sample, background signal subtracted.

**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