

Product datasheet

# Alcohol dehydrogenase Assay Kit ab272518

1 Image

Overview

<b>Product name</b>	Alcohol dehydrogenase Assay Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell Lysate, Tissue Lysate
<b>Assay type</b>	Quantitative
<b>Range</b>	0.4 mU/ml - 80 mU/ml
<b>Product overview</b>	<p>Alcohol Dehydrogenase Assay Kit (ab272518) is a non-radioactive, colorimetric ADH assay is based on the reduction of the tetrazolium salt MTT in a NADH coupled enzymatic reaction to a reduced form of MTT which exhibits an absorption maximum at 565 nm. The increase in absorbance at 565 nm is directly proportional to the enzyme activity.</p> <p><b>Fast and Sensitive:</b> Linear detection range (20 µL sample): 0.4 to 80 U/L for 30 min reaction. Detection Limit of 0.1 U/L for 120 minute reaction.</p> <p><b>Linear detection range (20 µL sample):</b> 0.4 to 80 U/L for 30 min reaction. Detection Limit of 0.1 U/L for 120 minute reaction.</p>
<b>Tested applications</b>	<b>Suitable for:</b> Functional Studies
<b>Platform</b>	Microplate

Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer	1 x 10ml
Calibrator	1 x 1.5ml
Diaphorase	1 x 120µl
MTT Solution	1 x 1.5ml
NAD Solution	1 x 1ml

<b>Components</b>	<b>100 tests</b>
Substrate (10% Ethanol)	1 x 1ml

**Relevance**

Alcohol dehydrogenase (Alcohol DH, ADH) (EC 1.1.1.1) is a group of seven dehydrogenase enzymes that occur in many organisms and facilitate the interconversion between alcohols and aldehydes or ketones with the reduction of NAD<sup>+</sup> to NADH. In humans and many other animals, they serve to break down alcohols which could otherwise be toxic; in yeast and many bacteria, some alcohol dehydrogenases catalyze the opposite reaction as part of fermentation.

**Cellular localization**

Cytoplasmic

**Applications**

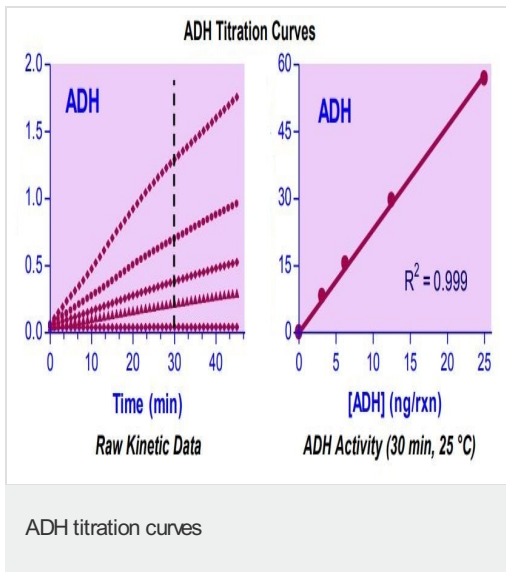
**The Abpromise guarantee**

Our [Abpromise guarantee](#) covers the use of ab272518 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
<b>Functional Studies</b>		Use at an assay dependent concentration.

**Images**



**Typical titration curves – data provided for demonstration purposes only.**

Example of Alcohol Dehydrogenase titration curves.

OD<sub>565</sub> readings (left) and calculated activity (right).

**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.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

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