

Alcohol Dehydrogenase Assay Kit ab102533

★★★★★ [1 Abreviews](#) [15 References](#) [2 Images](#)

Overview

Product name	Alcohol Dehydrogenase Assay Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Cell culture media
Assay type	Enzyme activity
Sensitivity	= 0.01 mU/well
Assay time	0h 40m
Product overview	<p>Abcam's Alcohol Dehydrogenase Assay Kit provides a convenient tool for sensitive detection of the Alcohol DH in a variety of samples. In the assay Alcohol DH will utilize isopropanol as a substrate leading to a proportional color development. The activity of ADH can be easily quantified colorimetrically ($\lambda = 450 \text{ nm}$). This assay detects ADH activity as low as 0.01 mU in samples.</p> <p>Visit our FAQs page for tips and troubleshooting.</p> <p>Alcohol dehydrogenase assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix - incubate for 3 min - analyze with microplate reader, incubate for 30 min to 2 hrs and analyze again
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K787 Alcohol Dehydrogenase Activity Colorimetric Assay Kit. K787-100 is the same size as the 100 test size of ab102533.</p> <p>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⁺ 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.</p>
Platform	Microplate reader

Properties

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

Components	100 tests
ADH Positive Control	1 vial
Assay Buffer XXXIV	1 x 25ml
Developer VIII	1 vial
Substrate XII	1 x 1ml
NADH Standard I	1 vial

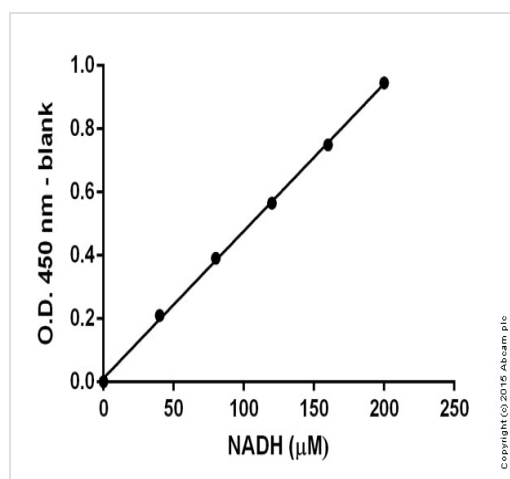
Relevance

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⁺ 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

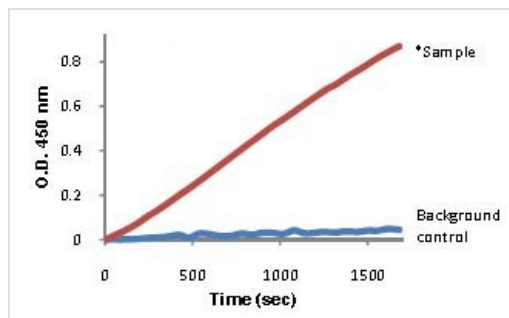
Cytoplasmic

Images



Colorimetric standard curve: mean of duplicates (+/-SD) with background readings subtracted.

Functional Studies - Alcohol Dehydrogenase
Detection Kit (ab102533)



Sample: Bovine Liver extraction (36µg protein)

Functional Studies - Alcohol Dehydrogenase
Detection Kit (ab102533)

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