

Alcohol Dehydrogenase Activity Assay

A plate based colorimetric assay to quantifying Alcohol Dehydrogenase (ADH)

Alcohol dehydrogenase plays a small metabolic role in humans but is a much more important contributor to the health and survivability of microorganisms such as yeast and bacteria. In humans ADH resides primarily in the liver where it detoxifies ethanol. Its role in microorganisms is more interesting as it is involved in anaerobic energy production, adaptability to toxic environments bioremediation. ADH generates NADH in the process of oxidizing alcohol, an important intermediate in energy production. AkrisBio's Alcohol Dehydrogenase Assay provides a simple, sensitive means of quantifying ADH activity in a variety of sample types with a lower limit of detection in the low microunit range.

Ordering Information

[Order Online »](#)

AKR-MA-0162	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com