

Ethanol Assay

A plate based colorimetric/fluorometric assay to measure ethanol concentration in a variety of samples.

Ethanol plays a role in various metabolic pathways within the human body. When consumed, it is metabolized primarily in the liver via oxidative metabolism, involving alcohol dehydrogenase and acetaldehyde dehydrogenase, converting ethanol through toxic acetaldehyde, to acetate and eventually CO₂ and water. Ethanol metabolism affects energy balance and nutrient utilization. The metabolism of ethanol takes priority over other macronutrients, leading to storage of excess carbohydrates and fats. Chronic alcohol consumption disrupts metabolic pathways, impacting liver function, lipid metabolism, and glucose regulation. AkrivisBio's Ethanol Assay provides a simple, sensitive method for precise measurement of ethanol in a variety of samples. In the assay, alcohol oxidase oxidizes ethanol forming H₂O₂ which is used to form resorufin with intense color and fluorescence with sensitivity to 0.05 nmol.

Ordering Information

[Order Online »](#)

AKR-MA-0131	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