

Alanine Aminotransferase (ALT) Assay

A plate based colorimetric/fluorometric assay to quantify Alanine aminotransferase activity in a variety of biological samples with a detection limit below 10 mU per well.

Alanine aminotransferase (ALT) is an enzyme also known as SGPT (serum glutamic pyruvic transaminase), and is found in serum and several tissues, predominantly in liver cells. High levels of ALT in the blood is an indication of liver damage or disease and is as part of a differential diagnostic test.

AkrivisBio's ALT Assay is a simple and sensitive assay to measure ALT enzyme activity in a variety of biological samples. In this assay, alanine and α -ketoglutarate are converted to pyruvate and glutamate in the presence of ALT. Pyruvate is then oxidized forming hydrogen peroxide, which is utilized by the peroxidase to oxidize ADHP thereby resulting in the formation of intense color or fluorescence measured at OD 570 nm or at Ex/Em = 535/587 nm respectively.

Ordering Information

[Order Online »](#)

AKR-MA-0116	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