

Aspartate Aminotransferase Activity Assay

A plate based colorimetric to measure AST activity in a variety of biological samples

Aspartate aminotransferase (AST), also known as serum glutamic oxaloacetic transaminase (SGOT), is an enzyme found in various tissues, with higher concentrations in the liver and heart. Measurement of AST levels is used to assess liver damage or disease. AkrivisBio's AST Activity Assay is a simple, sensitive way to measure AST activity in a wide range of biological samples with a detection limit of less than 0.02 mU per well. In this assay, AST catalyzes the conversion of aspartate and α -ketoglutarate to produce oxaloacetate and glutamate. Glutamate is then oxidized by glutamate dehydrogenase forming NADH in the process. NADH reduces WST, a readily reducible nearly colorless tetrazolium to a highly colored formazan measured at OD 450 nm.

Ordering Information

[Order Online »](#)

AKR-MA-0120	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