

Glutamate Assay

A plate based colorimetric assay to measure the concentration of glutamate in samples as low as 0.2 nmoles per well.

Glutamate, a non-essential amino acid, plays several important metabolic roles in the body. As a neurotransmitter, it facilitates communication between nerve cells, contributing to brain function and cognition. Glutamate is also a key molecule in protein synthesis, aiding in the production of other amino acids. It also serves as a precursor for glutathione, a potent antioxidant that helps protect cells from oxidative damage. AkrivisBio's Glutamate assay is a simple, sensitive means of assessing glutamate content in a variety of tissues with an ability to detect less than 0.2 nmoles per well. In the assay, Glutamate is oxidatively converted to α -ketoglutarate by glutamate dehydrogenase with the conversion of NAD to NADH. NADH then reduces a nearly colorless tetrazolium to a highly colored formazan measured at OD 450 nm. Glutamate assay is useful for studying neurotransmission, assessing neurological disorders, and evaluating glutamate levels in food products to monitor quality or detect contamination.

Ordering Information

[Order Online »](#)

AKR-MA-0133	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