

Glutamate Assay Kit

Glutamate is a non-essential amino acid that serves as an important neurotransmitter in the mammalian brain and has a key role in cellular metabolism. Excess glutamate levels in the brain can cause cell injury and death, leading to neurological diseases. Our Glutamate Assay Kit is a quantitative, fluorometric assay that uses glutamate specific enzymes to generate hydrogen peroxide. An ADHP probe is oxidized by hydrogen peroxide to generate fluorescent Resorufin, which correlates to the level of glutamate in the sample. Glutamate levels in an unknown sample are calculated based on a glutamate standard curve.

Ordering Information

[Order Online »](#)

CBI-STA-674	200 assays
-------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Detection sensitivity limit of 0.3 μM glutamate
- Suitable for use with lysates, cell culture supernatants, serum, plasma, and urine

Handling & Storage

Use/Stability Upon receipt, store the 10X Assay Buffer at 4°C. Aliquot and store all other components at -20°C. Avoid multiple freeze/thaw cycles. ADHP is light sensitive and must be stored accordingly.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescent detection

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](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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