

Glutamine Assay

A plate based colorimetric assay to quantify glutamine from a variety of biological samples with a sensitivity in the low nanomole range.

Glutamine, a non-essential amino acid, plays several important metabolic roles. Glutamine demand often outstrips production capacity and must be supplied by the diet. While important for protein synthesis, other roles are equally important. Glutamine is readily converted to glucose, particularly when intense physical activity or other stress increases glucose demand. Glutamine is both a source of nitrogen for synthesis of other molecular classes such as nucleotides and as the primary pathway for ammonia detoxification where it is used by the liver to produce urea allowing for excretion of excess amine/ammonia. Glutamine is essential for immune cells, particularly lymphocytes and macrophages, bolstering the immune response. Finally, glutamine is important to the gastrointestinal tract, providing the primary source of energy for the rapidly dividing cells of the intestinal lining. AkrivisBio's Glutamine Assay provides a simple sensitive means of quantifying glutamine from a variety of biological samples with a sensitivity in the low nanomole range.

Ordering Information

[Order Online »](#)

AKR-MA-0164	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