

# L-Asparaginase Activity Assay

A plate-based colorimetric/fluorometric assay to measure asparaginase activity in various biological samples.

Asparaginase catalyzes the hydrolysis of L-asparagine to L-aspartate and ammonia. Asparaginase does not occur naturally in humans but is found in bacteria, plants and some animals. Since asparagine is an essential amino acid in certain cell types, including some cancers, controlling the circulating levels of asparagine is a way of inhibiting protein synthesis needed for rapid growth. AkrisBio's Asparaginase Activity Assay provides a simple, sensitive means of measuring asparaginase activity in biological samples.

## Ordering Information

[Order Online »](#)

AKR-MA-0159	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](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)