

Trypsin Activity Assay

A plate based colorimetric assay to quantify Trypsin activity in a wide range of samples.

Trypsin is a serine protease that plays a crucial role in protein digestion and various other biological processes. It is produced in the pancreas in an inactive proenzyme form and then secreted into the small intestine where the proenzyme is activated by enterokinase. AkrivisBio's Trypsin Assay is a kinetic assay for the quantification of trypsin activity in a wide range of samples. In the assay, trypsin cleaves a trypsin substrate (GPK-pNA) to generate p-nitroaniline, which is detected at OD 405 nm. The intensity of the color is proportional to the p-nitroaniline content, trypsin activity. The kit can detect trypsin activity as low as 0.05 - 0.1 mU per sample.

Ordering Information

[Order Online »](#)

AKR-MA-0128	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°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