

α -Amylase Activity Assay

A plate based colorimetric assay to measure the enzymatic activity of α -amylase in samples.

Amylases hydrolyze starch and glycogen into maltose and oligosaccharides. α -Amylase is the major form of amylase that is found in humans. It is a calcium metallo-enzyme and is found in various tissues and body fluids including saliva, pancreatic juice, and blood. Its activity is commonly measured to assess the digestive function and diagnose various pathological conditions such as salivary gland trauma or inflammation, mumps, pancreatitis and renal failure. AkrivisBio's α -Amylase Activity Assay is a simple, sensitive means of measuring amylase activity in various sample types. In the assay, amylase cleaves the artificial substrate, ethylidene-pNP-G7 to small oligosaccharides. The oligosaccharides are subsequently hydrolyzed by α -glucosidase to release nitrophenol, which is measured at absorbance 405 nm. The kit can detect amylase activity at levels below 1 mU.

Ordering Information

[Order Online »](#)

AKR-MA-0138	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