

α -Galactosidase Activity Assay

A plate based fluorometric assay to measure the enzymatic activity of α -galactosidase in biological samples.

α -Galactosidase hydrolyzes α -galactosyl residues found in glycolipids, polysaccharides and glycoproteins. In mammals, α -Galactosidase is produced by the pancreas and released into the small intestine, where it hydrolyzes galactosides from dietary sources. AkrivisBio's α -Galactosidase Activity Assay is a simple, sensitive means of monitoring α -Galactosidase activity in biological samples. In the assay a synthetic substrate is acted on by α -Galactosidase releasing a fluorophore measured at Ex/Em = 360/445 nm. The α -galactosidase activity assay is a valuable research tool to study genetic disorders involving α -galactosidase deficiencies such as Fabry disease, and the development of potential therapeutic agents that modulate α -galactosidase activity.

Ordering Information

[Order Online »](#)

AKR-MA-0142	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 Fluorescent 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