

Lipase Activity Assay

A plate based colorimetric/fluorometric assay to measure lipase activity in a variety of biological samples.

Lipases play a pivotal role in the digestion, absorption, and metabolism of dietary fats in organisms. In the human digestive system, lipases break down triglycerides into fatty acids and glycerol, enabling efficient nutrient absorption in the intestine. Lipase is a key regulator of energy balance. By releasing fatty acids from adipose tissue during periods of energy demand, lipases enable the body to access stored energy reserves. This is particularly crucial during fasting or physical activity when glucose availability is limited. Lipase activity influences lipid-related disorders such as obesity, diabetes, and cardiovascular diseases. Dysregulation of lipase function can lead to abnormal lipid profiles and contribute to the development of these conditions. AkrivisBio's Lipase Activity Assay I is a simple sensitive way to measure lipase activity in a variety of biological samples.

Ordering Information

[Order Online »](#)

AKR-MA-0149	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