

Sulfatase Activity Assay

A plate based colorimetric assay to determine sulfatase activity in a variety of biological samples with a sensitivity below 1 mU per well.

Sulfatases are a group of 17 enzymes which catalyze the removal of the sulfate group from a variety of steroids and complex carbohydrates such as glycosaminoglycans. Deficiencies in sulfatase activity leads to lysosomal storage diseases where the sulfated molecules accumulate, causing functional impairment and tissue damage. The sulfatases also play a role in the regulation of hormone activity, deconjugating the sulfated forms of hormones like estrogen and DHEA which influences bioavailability and receptor binding. AkrivisBio's Sulfatase Activity Assay is a simple, sensitive means of assessing sulfatase activity in a variety of biological samples with a sensitivity below 1 mU per well.

Ordering Information

[Order Online »](#)

AKR-MA-0163	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