

Acetate Assay

A plate based colorimetric assay to quantify acetate levels in various samples.

Acetate, a molecule derived from various metabolic pathways, plays an important role in various biological processes, including energy metabolism and the synthesis of lipids. Additionally, acetate is a signaling molecule regulating gene expression and cellular processes. In AkrivisBio's Acetate Assay, acetate is phosphorylated, converting ATP to ADP. ADP is used to convert glucose to glucose-6-phosphate, which in turn is used to convert NAD to NADH. The NADH then reduces a nearly colorless tetrazolium to a highly colored formazan measured at OD 450 nm. The level of acetate quantitation is <1 nmole/well, equivalent to concentrations in the order of 20 μ M in tested samples. Acetate assay helps in understanding acetate levels assessing microbial activity, and evaluating metabolic changes in various physiological and pathological conditions.

Ordering Information

[Order Online »](#)

AKR-MA-0130	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