

Uric Acid Assay

A plate based colorimetric/fluorometric assay to measure uric acid levels in a variety of biological samples with a sensitivity below 1 nmole.

Uric acid, a byproduct of purine metabolism, has both positive and negative aspects in metabolism. While elevated levels are associated with gout, a painful joint condition, uric acid serves essential functions, acting as an antioxidant, neutralizing harmful free radicals, and protecting cells from oxidative stress. It contributes to the regulation of blood pressure by modulating the production of nitric oxide, a vasodilator. Despite its physiological benefits, excessively high uric acid levels can lead to gout and other metabolic disorders such as insulin resistance, obesity, and cardiovascular disease. AkrisBio's Uric Acid assay is a simple, sensitive means of measuring uric acid levels in a variety of biological samples with a sensitivity below 1 nmole.

Ordering Information

[Order Online »](#)

AKR-MA-0151	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