

Creatinine Assay

A plate based colorimetric/fluorometric assay to measure creatinine concentration in a variety of biological samples. The assay can quantify less than 1 nmole per well by absorbance and below 0.1 nmole per well by fluorescence.

Creatinine is a breakdown product of muscle and protein metabolism. Creatinine is produced and excreted at a constant rate, and as such is used as an internal control for the rate of formation or breakdown of a variety of metabolic intermediates. Since urine production rates can vary due to a number of factors, the creatinine concentration corrects for this and is used to normalize urine concentrations of other compounds. Creatinine is excreted via glomerular filtration and secretion from the proximal tubules. Blood creatinine levels increase if glomerular filtration decreases and are used to determine the estimated GFR (eGFR). AkriBio's Creatinine Assay Kit provides a simple, sensitive measure of creatinine concentration in a variety of biological samples. In the assay, creatinine broken down with the eventual formation of hydrogen peroxide which is used by peroxidase to convert ADHP to the highly colored and fluorescent resorufin. The assay can quantify less than 1 nmole per well by absorbance and below 0.1 nmole per well by fluorescence.

Ordering Information

[Order Online »](#)

AKR-MA-0121	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