

D-Lactate Assay

A plate based colorimetric assay to measure D-lactate levels in a variety of biological samples.

D-lactate, an enantiomer of the common L-lactate, is primarily produced by gut bacteria. When the balance of gut bacteria is disrupted, as seen in conditions such as short bowel syndrome, intestinal bacterial overgrowth, or certain surgical interventions, D-lactate levels can rise abnormally. This elevation can lead to D-lactic acidosis, a condition characterized by an accumulation of D-lactate in the blood, which can lead to neurological symptoms. D-lactate measurements can be used as an indicator of intestinal health and bacterial overgrowth, especially in patients with altered gut anatomy or function. AkrisBio's D-Lactate Assay is a simple, sensitive way to measure D-Lactate levels in a variety of biological samples.

Ordering Information

[Order Online »](#)

AKR-MA-0148	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