

L-Carnitine Assay

A plate based colorimetric/fluorometric assay to measure free L-Carnitine in biological samples.

Carnitine is a critical transport molecule which shuttles fatty acids into the mitochondria where they are utilized to make acetyl CoA, a key part of the TCA cycle. In the reverse direction carnitine shuttles toxic by-products of fatty acid metabolism out of the mitochondria, minimizing oxidative stress and oxidative damage that would otherwise occur. L-Carnitine is synthesized from methionine and lysine. Beyond its highly important shuttle function, it also enhances insulin sensitivity and cellular glucose uptake. AkrivisBio's L-Carnitine Assay is a simple, sensitive means of measuring free L-Carnitine in biological samples.

Ordering Information

[Order Online »](#)

AKR-MA-0161	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