

Succinate Assay

A plate based colorimetric assay to quantify succinate at levels below 25 μM in a variety of sample types.

Succinate is an important metabolite in the tricarboxylic acid cycle. Succinate has also emerged as a signaling molecule in various cellular processes, including inflammation, hypoxia responses, and regulation of gene expression. In pathological conditions such as ischemia, it accumulates, impairing mitochondrial function and promoting oxidative stress and tissue damage. AkrivisBio's Succinate assay is a simple, sensitive method for the accurate determination of succinate in a variety of samples at levels below 25 μM . In the assay, succinate is converted to succinyl CoA with the formation of an intermediate, which then undergoes a series of reactions to generate NADH. The NADH formed then reduces nearly colorless tetrazolium to a highly colored formazan measured at OD 450 nm.

Ordering Information

[Order Online »](#)

AKR-MA-0134	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