

Calcium Assay

A plate-based colorimetric assay to measure 0-1 µg of calcium levels in samples.

Calcium is one of the most important metal ions that is essential in the regulation of diverse cellular processes in all living organisms. They function as a key second messenger through its sequestration and release into the cytoplasm via calcium channels thereby regulating various cell signaling pathways. AkrivisBio's Calcium Assay provides a simple and direct method for measuring Ca²⁺ concentration in a variety of samples. In this assay, Ca²⁺ binds to alkaline O-Cresolphthalein Complexone to form a chromogen measured at OD 575 nm. The intensity of the color is directly proportional to the calcium concentration in the sample.

Ordering Information

[Order Online »](#)

AKR-MA-0104	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°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