

NAD/NADH Assay

A plate based colorimetric assay to measure NAD⁺ and NADH, and their ratio in various samples.

The NAD/NADH redox couple is essential for maintaining cellular redox homeostasis. NADH is the primary vehicle whereby reducing equivalents are converted to ATP in the cell. With thousands of oxidoreductases in the cell, each with its own set of optimal conditions, the assay of nicotinamide nucleotides becomes a useful very broad index of the overall health of the cell. AkrivisBio's NAD/NADH Assay provides a simple, sensitive means of detecting NAD and NADH. The ability to measure these concentrations allows the determination of the NAD/ NADH ratio, which gives important insight into the overall redox state of the cell. The assay is specific for the NAD/NADH couple and not NADP/ NADPH, which are used primarily for unrelated anabolic reactions.

Ordering Information

[Order Online »](#)

AKR-MA-0114	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