

NADP-NADPH Assay

A plate based colorimetric assay to measure NADP and/or NADPH levels in a wide range of biological samples.

NADP and NADPH play critical roles in metabolism and redox reactions. NADP acts as a coenzyme in various enzymatic reactions, particularly photosynthesis and oxidative stress defense. NADPH, the reduced form of NADP, is essential for biosynthetic reactions, such as fatty acid and nucleotide synthesis. Moreover, NADPH serves as a potent reducing agent, protecting cells from oxidative damage and maintaining redox homeostasis. The dynamic interplay between NADP and NADPH is indispensable for cellular function and overall metabolic balance. AkrivisBio's NADP/NADPH Assay is a simple, sensitive colorimetric method for measuring NADP and/or NADPH as low as 100 picomole levels in a wide range of biological samples. In this assay, NADP in the sample is used by an NADP specific alcohol dehydrogenase to form NADPH, which then reduces a nearly colorless tetrazolium salt to a highly colored formazan measured at 450nm. The assay does not recognize NAD⁺/NADH.

Ordering Information

[Order Online »](#)

AKR-MA-0126	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