

NAD⁺/NADH Assay Kit (Fluorometric)

Our NAD⁺/NADH Assay Kit detects NAD⁺ and NADH in cell and tissue lysates. Total NAD⁺/NADH can be detected or samples can be treated with an acid or base treatment to specifically detect NAD⁺ or NADH. This assay uses an enzymatic cycling reaction that reduces NAD⁺ to NADH, which then reacts with a fluorometric probe and is detected with a fluorescence plate reader at 530-570nm excitation and 590-600nm emission. NAD⁺/NADH levels in unknown samples are calculated based on the provided NAD⁺ standard curve.

Ordering Information

[Order Online »](#)

CBI-MET-5030	100 assays
--------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Detects NAD⁺, NADH, or total NAD⁺/NADH
- Compatible with cell or tissue lysates
- NAD⁺ standard included

Handling & Storage

Use/Stability Upon receipt, store the NAD Cycling Enzyme and NAD⁺ Standard at -80°C. Store the Fluorometric Probe and Electron Mediator at -20°C. Store the remaining components at 4°C.

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescent detection

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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)