

ϵ -NADP⁺

ϵ -NADP⁺ is a fluorescent analogue of β -NADP⁺ with λ_{exc} 300 nm and λ_{em} 410 nm. It can replace β -NADP⁺ as a substrate for corresponding enzymes and is thus useful in examining the binding and coenzyme activity of β -NADP⁺. The potential metabolite ϵ -ADPRP (Prod. No. BLG-E016) is available as well.

*BLG-N024-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

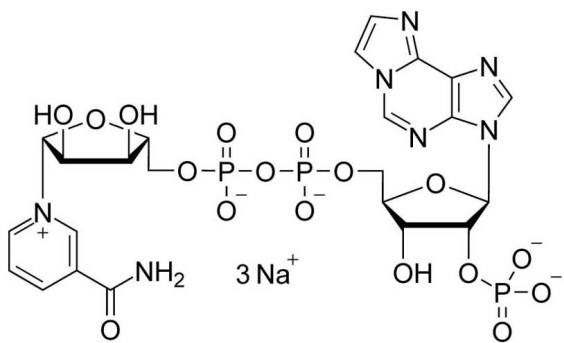
Ordering Information

[Order Online »](#)

BLG-N024-05	5 μ mol
BLG-N024-25	5x5 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling Protect from light.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name β -Nicotinamide-1,N6-ethenoadenine dinucleotide phosphate . sodium salt

CAS 104809-28-1

Couple Type Modified nucleotides

Formula $C_{23}H_{28}N_7O_{17}P_3$ (free acid)

MW 767.4 (free acid)

Purity > 95% HPLC

Quantity 5 μ mol = ~3.8 mg

Solubility Soluble in water (\geq 27mM).

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