

ϵ -NAADP⁺

ϵ -NAADP⁺ is a fluorescent analogue of the Ca²⁺-mobilizing agent nicotinic acid adenine dinucleotide phosphate (β -NAADP, Prod. No. [BLG-N018](#)) in which both the N¹ and the N⁶ nitrogen atoms in the adenine nucleobase are connected by an etheno bridge forming a tricyclic ring system.

ϵ -NAADP is a fluorescent analogue of the Ca²⁺-releasing messenger β -NAADP with λ_{exc} 275 nm and λ_{em} 410 nm. ϵ -NAADP activates Ca²⁺ release from sea urchin egg homogenates with a half-maximal effective concentration of 5 μ M (Lee & Aarhus 1998).

BLG-N019-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

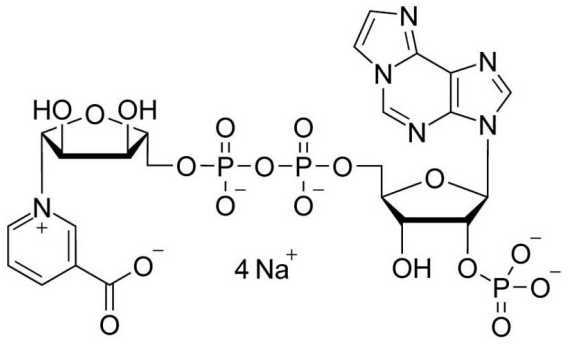
Ordering Information

[Order Online »](#)

BLG-N019-05	5 μ mol
BLG-N019-25	5x5 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Nicotinic acid-1,N6-ethenoadenine dinucleotide phosphate . tetrasodium salt

CAS 217313-72-9

Couple Type Modified nucleotides

Formula $C_{23}H_{27}N_6O_{18}P_3$ (free acid)

MW 768.4 (free acid)

Purity > 95% HPLC

Quantity 5 µmol = ~3.8 mg

Solubility Soluble in water (≥ 20 mM).

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