

β -NAADP⁺

β -NAADP⁺ is a potent stimulator of intracellular Ca²⁺ mobilization in a variety of cell types, with its mechanism of action being distinct from those of cyclic adenosine diphosphate ribose (cADPR, Prod. No. [BLG-C005](#)) and inositol-1,4,5-triphosphate (IP₃). The function of β -NAADP was first demonstrated in sea urchin egg homogenates.

BLG-N018-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

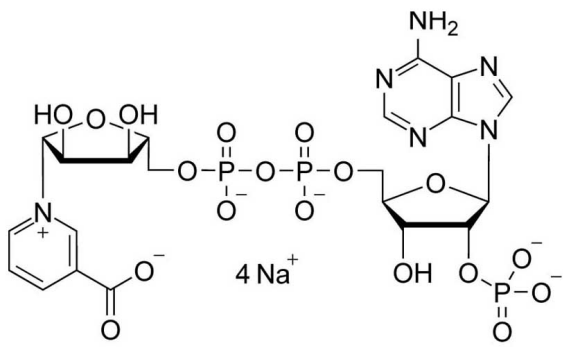
Ordering Information

[Order Online »](#)

| | |
|-------------|----------------|
| BLG-N018-10 | 10 μ mol |
| BLG-N018-50 | 5x10 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 adenine dinucleotide phosphate . tetrasodium salt

CAS 5502-96-5

Couple Type Modified nucleotides

Formula $C_{21}H_{27}N_6O_{18}P_3$ (free acid)

MW 744.4 (free acid)

Purity > 95% HPLC

Quantity 10µmol≈7.4mg

Solubility Soluble in water.

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