

6-AE-NAD⁺

functionalized for conjugation with markers
and dyes or for immobilization

6-AE-NAD⁺ is an analogue of β -NAD⁺ which is suitable as a ligand in affinity chromatography or for coupling of various labelling compounds including fluorophores.

*BLG-N013-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

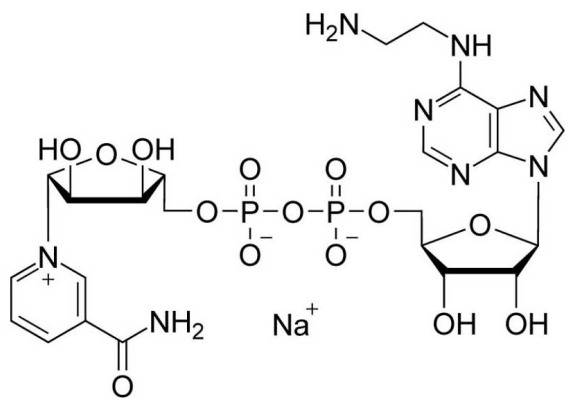
Ordering Information

[Order Online »](#)

BLG-N013-05	5 μ mol
BLG-N013-25	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 β -Nicotinamide-N6-(2-aminoethyl)adenine dinucleotide . sodium salt

CAS 59587-50-7

Couple Type Modified nucleotides

Formula $C_{23}H_{32}N_8O_{14}P_2$ (free acid)

MW 706.5 (free acid)

Purity > 95% HPLC

Quantity 5 μ mol \approx 3.5mg

Solubility Soluble in water (>58mM).

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