

8-BuT-NAD⁺

Analogue of β -NAD⁺, useful for screening of NAD receptor proteins.

*BLG-N036-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

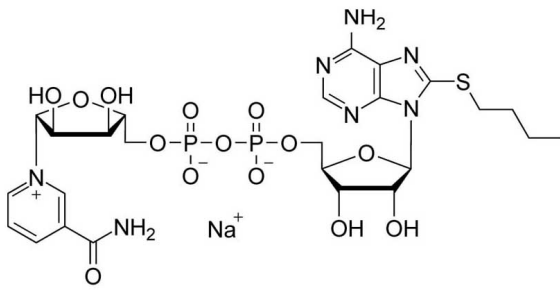
Ordering Information

[Order Online »](#)

BLG-N036-01	1 μ mol
BLG-N036-05	5x1 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-8-butylthioadenine dinucleotide . sodium salt

CAS 2022926-17-4

Couple Type Modified nucleotides

Formula $C_{25}H_{35}N_7O_{14}P_2S$ (free acid)

MW 751.6 (free acid)

Purity > 95% HPLC

Technical Info / Product Notes This product and/or its use with the asPARP technology are subject to patent application US2016299141 (A1) and foreign equivalents. Please visit www.biolog.de for further information.



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)