

# 8-Br-NAD<sup>+</sup>

8-Br-NAD<sup>+</sup> is an analogue of  $\beta$ -NAD<sup>+</sup> that can be utilized as a substrate by the ADP-ribosyl cyclase CD38 to produce the cADPR antagonist 8-Br-cADPR (Prod. No. [BLG-B065](#)).

*BLG-N017-50 (5 x 10  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

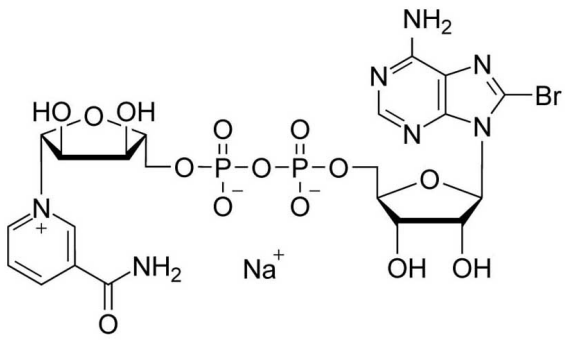
## Ordering Information

[Order Online »](#)

BLG-N017-50	5x10 $\mu$ 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**  $\beta$ -Nicotinamide-8-bromoadenine dinucleotide . sodium salt

**CAS** 2022926-16-3

**Couple Type** Modified nucleotides

**Formula**  $C_{21}H_{26}BrN_7O_{14}P_2$  (free acid)

**MW** 742.3 (free acid)

**Purity** > 95% HPLC

**Quantity** 10 $\mu$ mol $\approx$ 7.4mg

**Solubility** Soluble in water (>36mM).

**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](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)