

# 8-Br-7-CH-NAD<sup>+</sup>

8-Br-7-CH-NAD<sup>+</sup> is an analogue of  $\beta$ -NAD<sup>+</sup> in which the hydrogen in position 8 of the adenine nucleobase is replaced by bromine. In addition, the nitrogen atom in position 7 of the adenine imidazole ring is replaced by carbon and hydrogen, respectively. A potential substrate, competitive inhibitor or regulator of enzymes that interact with  $\beta$ -NAD<sup>+</sup>.

*BLG-N016-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

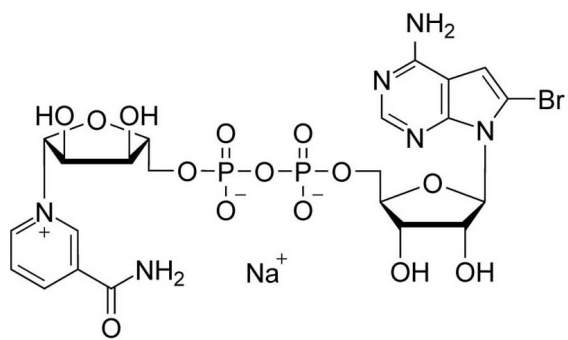
## Ordering Information

[Order Online »](#)

BLG-N016-01	1 $\mu$ mol
BLG-N016-05	5x1 $\mu$ 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-bromo-7-deazaadenine dinucleotide . sodium salt

**CAS** 189876-09-3

**Couple Type** Modified nucleotides

**Formula**  $C_{22}H_{27}BrN_6O_{14}P_2$  (free acid)

**MW** 741.3 (free acid)

**Purity** > 95% HPLC

**Quantity** 1  $\mu$ mol = ~0.7 mg

**Solubility** Soluble in water ( $\geq$  8 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](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)