

# carba-NAD<sup>+</sup>

**NAD<sup>+</sup> mimic with improved hydrolytic stability against adenosine diphosphate ribosyl transferase (ART) enzymes and NAD<sup>+</sup> glycohydrolases**

*BLG-C403-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

## Ordering Information

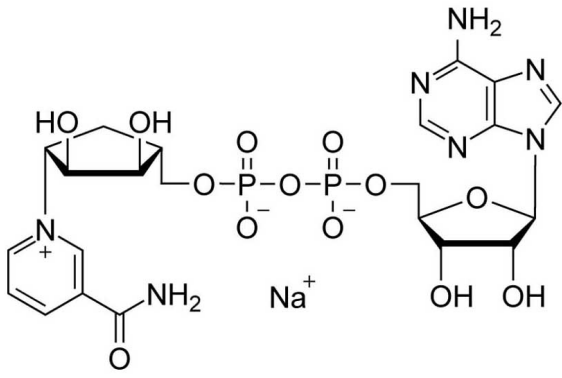
[Order Online »](#)

BLG-C403-01	1 $\mu$ mol
BLG-C403-05	5x1 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)

carba-NAD<sup>+</sup> is a NAD<sup>+</sup> mimic with improved hydrolytic stability against adenosine diphosphate ribosyl transferase (ART) enzymes and NAD<sup>+</sup> glycohydrolases. carba-NAD<sup>+</sup> was shown to inhibit activity of the cholera toxin A ART enzyme and NAD<sup>+</sup>-consuming enzymes Sirt3/5, but was ineffective against PARP1 and NAD<sup>+</sup> glycohydrolase. carba-NAD<sup>+</sup> is a substrate of NAD<sup>+</sup>-handling enzymes like alcohol dehydrogenases and is suitable as hydrolysis-resistant NAD<sup>+</sup>-mimic for protein crystallizations. Robust cofactor for enzyme-based glucose sensing with glucose dehydrogenases and fluorescence lifetime detection.



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Carba-β-nicotinamide adenine dinucleotide . sodium salt

**CAS** 112345-60-5

**Formula**  $C_{22}H_{29}N_7O_{13}P_2$  (free acid)

**MW** 661.5 (free acid)

**Purity** > 95% HPLC

**Quantity** 0.1 μmol / ~ 0.7 mg



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)