

carba-NAD⁺

NAD⁺ mimic with improved hydrolytic stability against adenosine diphosphate ribosyl transferase (ART) enzymes and NAD⁺ glycohydrolases

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

Ordering Information

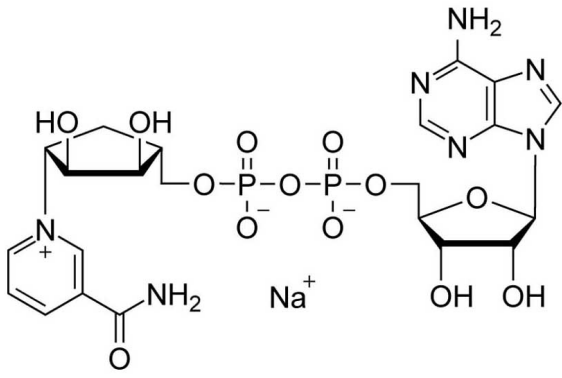
[Order Online »](#)

BLG-C403-01	1 μ mol
BLG-C403-05	5x1 μ mol

Manuals, SDS & CofA

[View Online »](#)

carba-NAD⁺ is a NAD⁺ mimic with improved hydrolytic stability against adenosine diphosphate ribosyl transferase (ART) enzymes and NAD⁺ glycohydrolases. carba-NAD⁺ was shown to inhibit activity of the cholera toxin A ART enzyme and NAD⁺-consuming enzymes Sirt3/5, but was ineffective against PARP1 and NAD⁺ glycohydrolase. carba-NAD⁺ is a substrate of NAD⁺-handling enzymes like alcohol dehydrogenases and is suitable as hydrolysis-resistant NAD⁺-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

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