

# BAD

Analogue of  $\beta$ -NAD<sup>+</sup> with improved stability

BLG-B301-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.

## Ordering Information

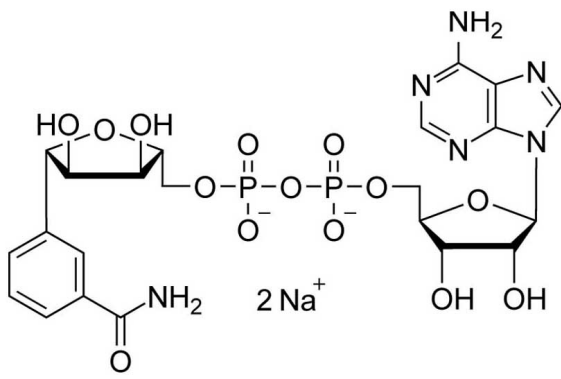
[Order Online »](#)

BLG-B301-01	1 $\mu$ mol
BLG-B301-05	5x1 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)

BAD is the biologically active metabolite of benzamide riboside and a potent inhibitor of inosine 5'-monophosphate dehydrogenase (IMPDH). It was also found to be a potent and competitive inhibitor of human NAD kinase ( $K_i = 90 \mu\text{M}$ ) and, at higher concentration and to a lower extent, of the bacterial enzyme. Furthermore, BAD is a non-hydrolysable substrate/inhibitor of NADases and PARP enzymes. BAD was used in co-crystallization experiments with NADases and PARP1.



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name  $\beta$ -Benzamide adenine dinucleotide . sodium salt

CAS 156724-91-3

Excitation Maximum 259 nm

Formula  $C_{22}H_{48}N_6O_{14}P_2$  (free acid)

MW 662.4 (free acid)

Purity > 95% HPLC

Quantity 1  $\mu$ mol / ~ 0.7 mg

**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)