

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

Analogue of  $\beta$ -NAD<sup>+</sup>, useful for screening of NAD receptor proteins.

*BLG-N037-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

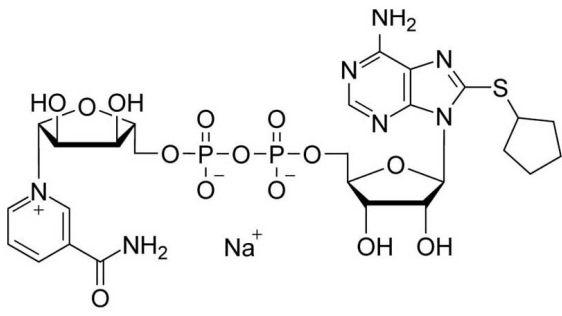
## Ordering Information

[Order Online »](#)

BLG-N037-01	1 $\mu$ mol
BLG-N037-05	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-cyclopentylthioadenine dinucleotide . sodium salt

**CAS** 2022926-10-7

**Couple Type** Modified nucleotides

**Formula**  $C_{26}H_{35}N_7O_{14}P_2S$  (free acid)

**MW** 763.6 (free acid)

**Purity** > 95% HPLC

**Technical Info / Product Notes** This product and/or its use with the asPARP technology are subject to patent application US2016299141 (A1) and foreign equivalents. Please visit [www.biolog.de](http://www.biolog.de) for further information.



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)