

# 6-Biotin-17-NAD<sup>+</sup>

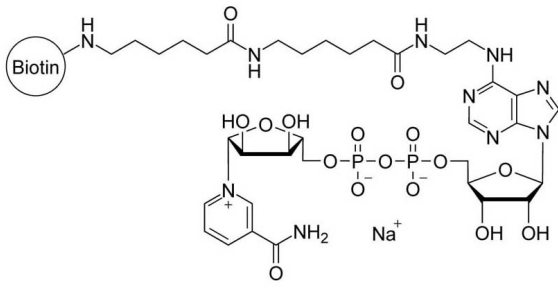
6-Biotin-17-NAD<sup>+</sup> is a biotinylated analogue of  $\beta$ -NAD<sup>+</sup> which can provide a non-radioactive alternative to radio-labelled  $\beta$ -NAD<sup>+</sup> for corresponding studies.

*BLG-N012-005 (5 x 0.1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

## Ordering Information [Order Online »](#)

BLG-N012-001	0.1 $\mu$ mol
BLG-N012-005	5x0.1 $\mu$ 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-N6-(2-(6-(6-[biotinyl]aminohexanoyl)aminohexanoyl)aminoethyl)adenine dinucleotide . sodium salt

**CAS** 1369508-04-2

**Couple Type** Modified nucleotides

**Formula**  $C_{45}H_{67}N_{12}O_{18}P_2S$  (free acid)

**MW** 1159.1 (free acid)

**Purity** > 95% HPLC

**Quantity** 0.1  $\mu$ mol $\approx$ 0.12mg

**Solubility** Soluble in water (>3.5mM).

**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)