

## Data Sheet

### **NAD<sup>+</sup>, Biotin-labeled**

Catalog #: 80610

Lot #: 220323

**Supplied as:** 250 µM aqueous solution**Stability:** Store solutions in aliquots at -80°C for up to one year.**References:**

1. Ying, W. (2008). *Antioxidants & Redox Signaling*. **10(2)**:179-206.
2. Ying, W. (2007). *Front Biosci*. **12**:1863-88.
3. Till, S., and Ladurner, A.G. (2009). *Front Biosci*. **14**:3246-3258.
4. Zhang, J. (1997). *Meth. Enzymol*. **280**: 255-65.

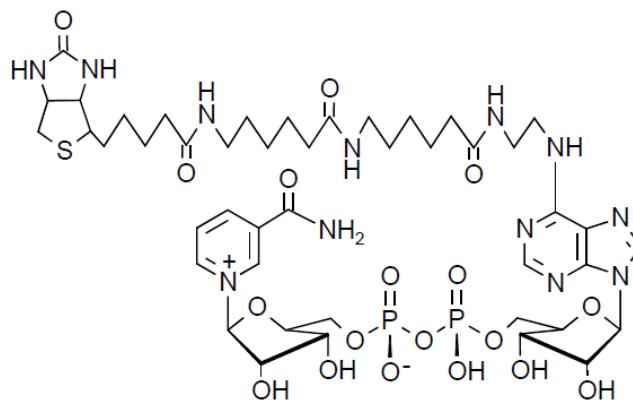
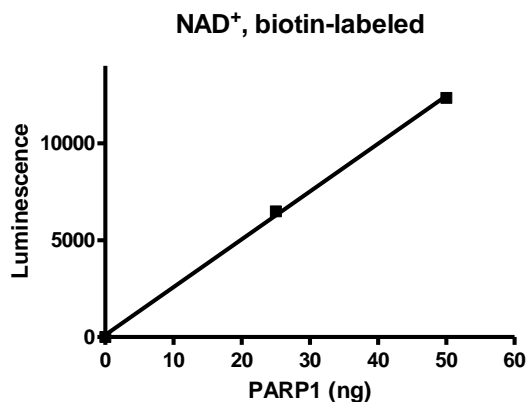
**Description:**

6-Biotin-17-NAD<sup>+</sup> (Nicotinamide adenine dinucleotide, or β-Nicotinamide- N6- (2- (6- [biotinyl] aminohexanoyl) aminoethyl) adenine dinucleotide). MW = 1158 (free acid).

**Purity:** ≥95% (HPLC)**Application:**

Useful for the study of PARP enzyme activity and DNA damage, or to purify ADP-ribosylated enzymes.

## **Quality Assurance**



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)