

# TREVIGEN® Product Data

*For Research Use Only. Not For Use In Diagnostic Procedures*

## 6-Biotin-17-NAD

**Catalog #:** 4670-500-01 **Size:** 500 µl

**Name:** 6-biotin-17-nicotinamide-adenine-dinucleotide

**Concentration:** 250 µM in water (Anion)

**Molecular Weight:** 1068 g/mole (Salt); 1045 g/mole (Anion)

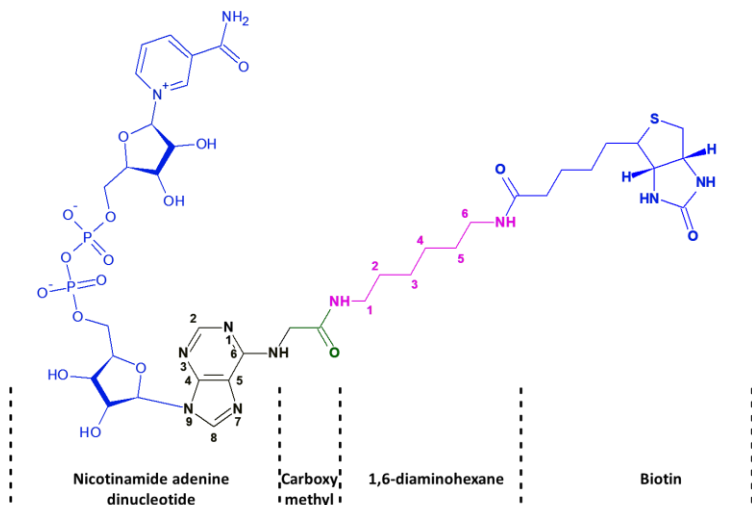
**Molecular Formula:** C<sub>39</sub>H<sub>56</sub>N<sub>11</sub>NaO<sub>17</sub>P<sub>2</sub>S (Salt); C<sub>39</sub>H<sub>56</sub>N<sub>11</sub>O<sub>17</sub>P<sub>2</sub>S (Anion)

**Extinction Coefficient:** 22,000 at 265 nm

**Purity:** Greater than 95% by HPLC

**Storage Conditions:** Store at -80°C. To avoid repeated freeze thaw cycles, prepare aliquots and store at -80°C.

**Structure:**



© 2016 Trevigen, Inc. All Rights Reserved. Trevigen is a registered trademark of Trevigen, Inc.

# TREVIGEN®

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: [info@trevigen.com](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

## Related Products Available From Trevigen

Catalog #	Description	Size
4676-096-K	Universal Chemiluminescent PARP Assay Kit/w Histone Coated Strip Wells	96 samples
4677-096-K	Universal Colorimetric PARP Assay Kit/w Histone Coated Strip Wells	96 samples

**6-Biotin-17-NAD**

Catalog #: 4670-500-01

Storage: -80 °C

**TREVIGEN®**  
1-800-873-8443