

# 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  
(C<sub>45</sub>H<sub>67</sub>N<sub>12</sub>O<sub>18</sub>P<sub>2</sub>SNa)

Concentration: 250 µM in water

Molecular Weight: 1180.08 g/mole (1158.09 g/mole as a free acid)

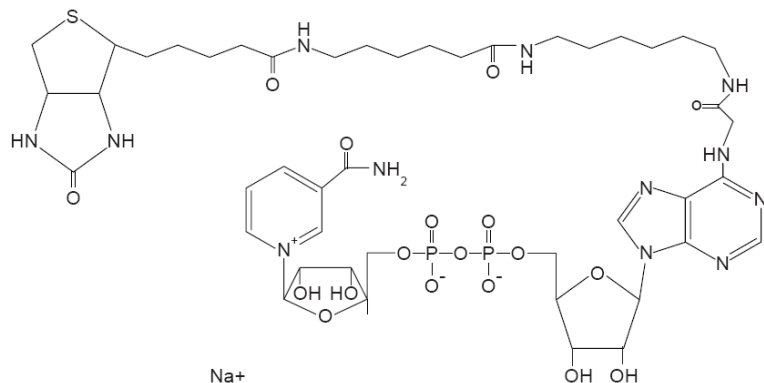
Extinction Coefficient: 22,000 at 265 nm

Purity: Greater than 95% by HPLC

Greater than 95% by <sup>31</sup>P NMR

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

### Structure:



**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)

**6-Biotin-17-NAD**  
Catalog #: 4670-500-01  
Storage: -80 °C  
**TREVIGEN®**  
1-800-873-8443