

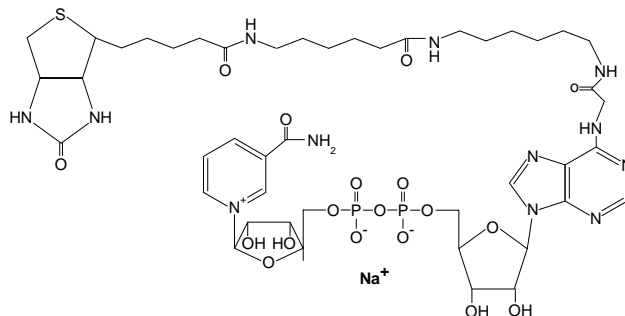
TREVIGEN™ Product Data

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 ₄₅ H ₆₇ N ₁₂ O ₁₈ P ₂ SNa)	
Concentration:	250 µM in water	
Molecular Weight:	1180.08 g/mole (1158.09 as free acid)	
Extinction coefficient:	22,000 at 265 nm	
Purity:	Greater than 95% by HPLC Greater than 95% by ³¹ P NMR	
Storage Conditions:	Store at -80°C. To avoid repeated freeze-thaw cycles, prepare aliquots and store at -80°C.	
Lot Specific Data:	See reverse	

Structure:



© 2000 Trevigen, Inc. All Rights Reserved. Trevigen is a trademarks of Trevigen, Inc.

v00530

TREVIGEN®

8405 Helgerman Court, Gaithersburg, MD 20877 USA
Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800
Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com