

D-Biotin

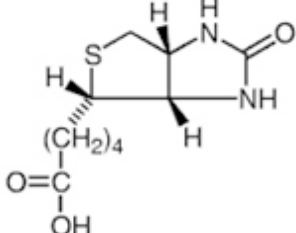
Catalog Number	Packaging Size
B001	1 g
B002	5 g

Storage upon receipt: -20°C

Introduction

Biotin (Known as Vitamin H or Coenzyme R) is a water-soluble B-complex vitamin (vitamin B7). D-biotin binds with high affinity to streptavidin and avidin. Biotin is a coenzyme for carboxylase enzymes, involved in the synthesis of fatty acids, isoleucine, and valine, and in gluconeogenesis.

Specifications

Molecular Formula:	C ₁₀ H ₁₆ N ₂ O ₃ S	
Molecular Weight:	244.31	
CAS Number:	58-85-5	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Biotinylation reagent

References:

1. [Easily reversible desthiobiotin binding to streptavidin, avidin, and other biotin-binding proteins: uses for protein labeling, detection, and isolation.](#)
Hirsch JD, Eslamizar L, Filanoski BJ, Malekzadeh N, Haugland RP, Beechem JM, Haugland RP
Anal Biochem (2002) 308:343-357
2. [A SPECTROPHOTOMETRIC ASSAY FOR AVIDIN AND BIOTIN BASED ON BINDING OF DYES BY AVIDIN.](#)
GREEN NM,
Biochem J (1965) 94:23C-24C