

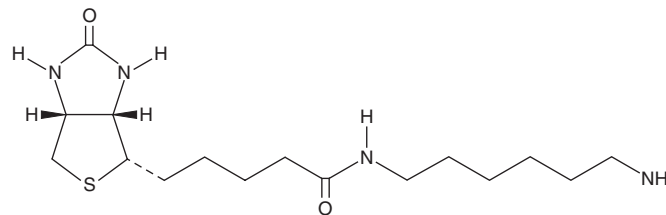
# PRODUCT INFORMATION



## Biotinyl hexylamine

Item No. 31226

**CAS Registry No.:** 65953-56-2  
**Formal Name:** (3a*S*,4*S*,6a*R*)-*N*-(6-aminohexyl)hexahydro-2-oxo-1*H*-thieno[3,4-*d*]imidazole-4-pentanamide  
**MF:** C<sub>16</sub>H<sub>30</sub>N<sub>4</sub>O<sub>2</sub>S  
**FW:** 342.5  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Biotinyl hexylamine is supplied as a solid. A stock solution may be made by dissolving the biotinyl hexylamine in the solvent of choice, which should be purged with an inert gas. Biotinyl hexylamine is slightly soluble in ethanol and chloroform.

### Description

Biotinyl hexylamine is a synthetic intermediate.<sup>1,2</sup> It has been used in the synthesis of avidin nucleic acid nano assemblies (ANANAS) and chemotherapeutic conjugates for the diagnosis of, and drug delivery to, tumors.

### References

1. Sabatino, G., Chinol, M., Paganelli, G., *et al.* A new biotin derivative-DOTA conjugate as a candidate for pretargeted diagnosis and therapy of tumors. *J. Med. Chem.* **46**(14), 3170-3173 (2003).
2. Bigini, P., Previdi, S., Casarin, E., *et al.* *In vivo* fate of avidin-nucleic acid nanoassemblies as multifunctional diagnostic tools. *ACS Nano* **8**(1), 175-187 (2014).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/18/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM