

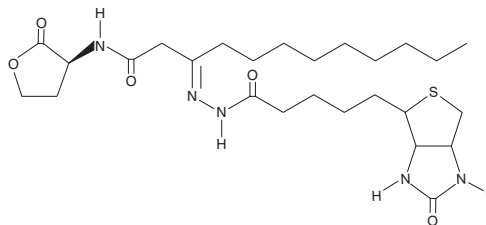
PRODUCT INFORMATION



N-dodecanoyl-L-Homoserine lactone-3-hydrazone-biotin

Item No. 9000710

Formal Name: N-[(3S)-tetrahydro-2-oxo-3-furanyl]-dodecanamide-hydrazone-biotin
Synonym: N-dodecanoyl-L-HSL-3-hydrazone-biotin
MF: $C_{26}H_{43}N_5O_5S$
FW: 537.7
Purity: $\geq 90\%$
UV/Vis.: λ_{max} : 238 nm
Supplied as: A solution in acetone
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

N-dodecanoyl-L-HSL-3-hydrazone-biotin is supplied as a solution in acetone. To change the solvent, simply evaporate the acetone under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of N-dodecanoyl-L-HSL-3-hydrazone-biotin in these solvents is approximately 3 and 5 mg/ml, respectively. While C₉-HSL is also soluble in ethanol and other primary alcohols, their use is not recommended as they have been shown to open the lactone ring.

Description

N-3-oxo-dodecanoyl-L-Homoserine lactone (Item No. 10007895) is a diffusible molecule involved in quorum sensing by bacteria. Homoserine lactones are also produced following proteolytic cleavage at a methionine residue by cyanogen bromide, as may be done during protein analysis. N-dodecanoyl-L-HSL-3-hydrazone-biotin is an affinity probe which allows the homoserine lactone molecule to be detected or captured through an interaction with an avidin-containing probe.

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, 05/07/2024

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