

PRODUCT INFORMATION



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

Item No. 9000711

Formal Name: N-[(3S)-tetrahydro-2-oxo-3-furanyl]-dodecanamide-hydrazone-fluorescein

Synonym: N-dodecanoyl-L-HSL-3-hydrazone-fluorescein

MF: C₃₇H₄₀N₄O₈S

FW: 700.8

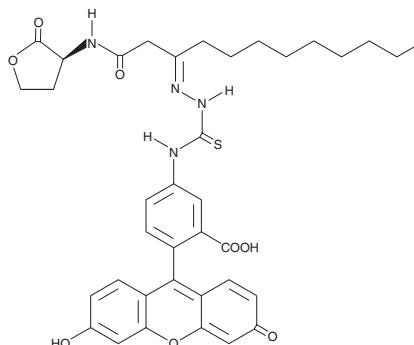
Purity: ≥95%

UV/Vis.: λ_{max}: 225, 277 nm

Supplied as: A crystalline 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

N-dodecanoyl-L-Homoserine lactone-3-hydrazone-fluorescein (N-dodecanoyl-L-HSL-3-hydrazone-fluorescein) is supplied as a crystalline solid. A stock solution may be made by dissolving the N-dodecanoyl-L-HSL-3-hydrazone-fluorescein in the solvent of choice, which should be purged with an inert gas. N-dodecanoyl-L-HSL-3-hydrazone-fluorescein is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml. While N-dodecanoyl-L-HSL-3-hydrazone-fluorescein 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-HSL (Item No. 10007895) is a diffusible molecule involved in quorum sensing by bacteria. HSLs 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-fluorescein is a fluorescently tagged probe which allows the HSL molecule to be detected using fluorescence-based techniques.

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, 04/23/2024

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