

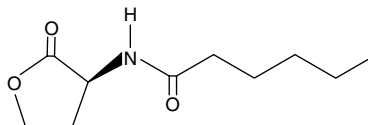
Product Information



N-hexanoyl-L-Homoserine lactone

Item No. 10007896

CAS Registry No.: 147852-83-3
Formal Name: N-[(3S)-tetrahydro-2-oxo-3-furanyl]-hexanamide
Synonym: C6-HSL
MF: C₁₀H₁₇NO₃
FW: 199.2
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that N-hexanoyl-L-homoserine lactone (C6-HSL) be stored as supplied at -20°C. It should be stable for at least two years.

C6-HSL is supplied as a crystalline solid. A stock solution may be made by dissolving the C6-HSL in organic solvents such as DMSO and dimethyl formamide. The solubility of C6-HSL in these solvents is approximately 30 mg/ml. While C6-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.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of C6-HSL can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of C6-HSL in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Quorum sensing is a regulatory system used by bacteria for controlling gene expression in response to increasing cell density. A promising field of study involves controlling bacterial infections by quenching their quorum sensing systems. The expression of specific target genes, such as transcriptional regulators belonging to the LuxIR family of proteins, is coordinated by synthesis of diffusible acylhomoserine lactone (AHL) molecules. C6-HSL is a small diffusible signaling molecule involved in quorum sensing, controlling gene expression, and affecting cellular metabolism.¹ The diverse applications of this molecule include regulation of virulence in general and in cystic fibrosis, infection prevention, slime and biofilm reduction in commercial agriculture and aquaculture industries, food spoilage prevention, and septicemia in fish.²⁻⁸

References

1. Gould, T.A., Herman, J., Krank, J., *et al. J. Bacteriol.* **188**(2), 773-783 (2006).
2. Morohoshi, T., Inaba, T., Kato, N., *et al. Journal of Bioscience and Bioengineering* **98**(4), 274-281 (2004).
3. Ulrich, R.L., Hines, H.B., Parthasarathy, N., *et al. J. Bacteriol.* **186**(13), 4350-4360 (2004).
4. Wang, L.-H., Weng, L.-X., Dong, Y.-H., *et al. J. Biol. Chem.* **279**(14), 13645-13651 (2004).
5. Yates, E.A., Philipp, B., Buckley, C., *et al. Infect. Immun.* **70**(10), 5635-5646 (2002).
6. Luo, Z.-Q., Su, S., and Farrand, S.K. *J. Bacteriol.* **186**(19), 5665-5672 (2003).
7. Chambers, C.E., Visser, M.B., Schwab, U., *et al. FEMS Microbiol. Lett.* **244**, 297-304 (2005).
8. Christensen, A.B., Riedel, K., Eberl, L., *et al. Microbiology* **149**, 471-483 (2003).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10007896

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 03/28/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com