

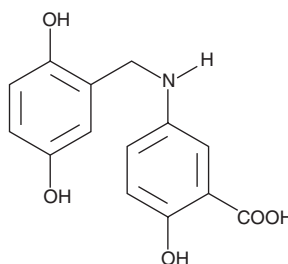
Product Information



Lavendustin C

Item No. 10010329

CAS Registry No.: 125697-93-0
Formal Name: 5-[[[(2,5-dihydroxyphenyl)methyl]amino]-2-hydroxy-benzoic acid
Synonyms: HDBA, NSC 666251
MF: C₁₄H₁₃NO₅
FW: 275.3
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that HDBA be stored as supplied at -20°C. It should be stable for at least two years.

HDBA is supplied as a crystalline solid. If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. HDBA is not stable and should be used immediately.

HDBA, a derivative of a *Streptomyces griseolavendus* butyl acetate extract, is a potent inhibitor of epidermal growth factor (EGF) receptor-associated tyrosine kinase with an IC₅₀ value of 0.012 μM.¹ HDBA also inhibits pp60^{c-src(+)} kinase and Ca²⁺ calmodulin-dependent kinase II with IC₅₀ values of 0.5 and 0.2 μM, respectively.² At a concentration of 10-150 μM, HDBA inhibits tyrosine kinase-associated neutrophil degranulation and superoxide generation.³

References

1. Onoda, T., Iinuma, H., Sasaki, Y., *et al.* Isolation of a novel tyrosine kinase inhibitor, lavendustin A, from *Streptomyces griseolavendus*. *J. Nat. Prod.* **52(6)**, 1252-1257 (1989).
2. O'Dell, T.J., Kandel, E.R., and Grant, S.G.N. Long-term potentiation in the hippocampus is blocked by tyrosine kinase inhibitors. *Nature* **353**, 558-560 (1991).
3. Burt, H.M., Jackson, J.K., and Salari, H. Inhibition of crystal-induced neutrophil activation by a protein tyrosine kinase inhibitor. *J. Leukoc. Biol.* **55**, 112-119 (1994).

Related Products

KN-62 - Item No. 13318 • Erbstatin Analog - Item No. 10010238 • Lavendustin A - Item No. 10010268 • RG-13022 - Item No. 10010309 • RG-14620 - Item No. 10010310

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

MATERIAL 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 Material 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, 03/21/2011

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