

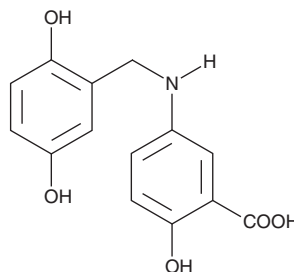
# PRODUCT INFORMATION



## 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:** C<sub>14</sub>H<sub>13</sub>NO<sub>5</sub>  
**FW:** 275.3  
**Purity:** ≥98%  
**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

Lavendustin C is supplied as a crystalline solid. A stock solution may be made by dissolving the lavendustin C in the solvent of choice, which should be purged with an inert gas. Lavendustin C is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of lavendustin C in these solvents is approximately 30 mg/ml.

### Description

Lavendustin C, a derivative of a *Streptomyces griseolavendus* butyl acetate extract, is a potent inhibitor of epidermal growth factor (EGF) receptor-associated tyrosine kinase with an IC<sub>50</sub> value of 0.012 μM.<sup>1</sup> Lavendustin C also inhibits pp60<sup>c-src(+)</sup> kinase and Ca<sup>2+</sup> calmodulin-dependent kinase II with IC<sub>50</sub> values of 0.5 and 0.2 μM, respectively.<sup>2</sup> At a concentration of 10-150 μM, lavendustin C inhibits tyrosine kinase-associated neutrophil degranulation and superoxide generation.<sup>3</sup>

### References

1. Onoda, T., Inuma, 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**(6344), 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**(1), 112-119 (1994).

#### 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, 01/26/2024

#### 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