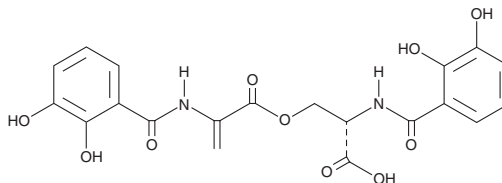


PRODUCT INFORMATION

N,N'-bis(2,3-Dihydroxybenzoyl)-O-L-seryl-L-dehydroalanine

Item No. 31779

CAS Registry No.: 73410-35-2
Formal Name: N-(2,3-dihydroxybenzoyl)-L-serine
 3-[2-[(2,3-dihydroxybenzoyl)amino]-2-propenoate]
MF: C₂₀H₁₈N₂O₁₀
FW: 446.4
Purity: ≥95%
Supplied as: A 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,N'-bis(2,3-Dihydroxybenzoyl)-O-L-seryl-L-dehydroalanine is supplied as a solid. A stock solution may be made by dissolving the N,N'-bis(2,3-dihydroxybenzoyl)-O-L-seryl-L-dehydroalanine in the solvent of choice, which should be purged with an inert gas. N,N'-bis(2,3-Dihydroxybenzoyl)-O-L-seryl-L-dehydroalanine is soluble in methanol and DMSO.

Description

N,N'-bis(2,3-Dihydroxybenzoyl)-O-L-seryl-L-dehydroalanine is a bis-catechol. It is a potential metabolite of the microbial metabolite N,N'-bis(2,3-dihydroxybenzoyl)-O-L-seryl-L-serine (Item No. 31778) based on the formation of dehydroalanine residues in serine-containing microbial peptides.¹ At the time N,N'-bis(2,3-dihydroxybenzoyl)-O-L-seryl-L-dehydroalanine was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available.

Reference

1. Arnison, P.G., Bibb, M.J., Bierbaum, G., *et al.* Ribosomally synthesized and post-translationally modified peptide natural products: Overview and recommendations for a universal nomenclature. *Nat. Prod. Rep.* **30**(1), 108-160 (2013).

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, 11/10/2022

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