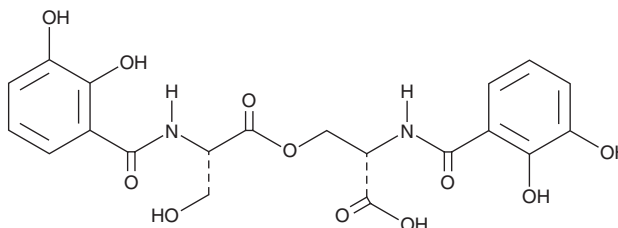


# PRODUCT INFORMATION

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

Item No. 31778

**CAS Registry No.:** 30414-15-4  
**Formal Name:** N-(2,3-dihydroxybenzoyl)-O-[N-(2,3-dihydroxybenzoyl)-L-seryl]-L-serine  
**MF:** C<sub>20</sub>H<sub>20</sub>N<sub>2</sub>O<sub>11</sub>  
**FW:** 464.4  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Bacterium/*Streptomyces* sp.



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-serine is supplied as a solid. A stock solution may be made by dissolving the N,N'-bis(2,3-dihydroxybenzoyl)-O-L-seryl-L-serine in the solvent of choice, which should be purged with an inert gas. N,N'-bis(2,3-Dihydroxybenzoyl)-O-L-seryl-L-serine is soluble in methanol and DMSO.

### Description

N,N'-bis(2,3-Dihydroxybenzoyl)-O-L-seryl-L-serine is a microbial metabolite and degradation product of the bacterial siderophore enterobactin that has been found in *Streptomyces* and has anticancer activity.<sup>1-3</sup> It inhibits invasion of 26-L5 murine colon cancer cells in a Matrigel™ assay with an IC<sub>50</sub> value of 2.7 μM.<sup>1</sup>

### References

1. Igarashi, Y., Iida, T., Fukuda, T., *et al.* Catechoserine, a new catecholate-type inhibitor of tumor cell invasion from *Streptomyces* sp. *J. Antibiot. (Tokyo)* **65**(4), 207-209 (2012).
2. O'Brien, I.G., Cox, G.B., and Gibson, F. Enterochelin hydrolysis and iron metabolism in *Escherichia coli*. *Biochim. Biophys. Acta* **237**(3), 537-549 (1971).
3. Raines, D.J., Moroz, O.V., Blagova, E.V., *et al.* Bacteria in an intense competition for iron: Key component of the *Campylobacter jejuni* iron uptake system scavenges enterobactin hydrolysis product. *Proc. Nat. Acad. Sci. USA* **113**(21), 5850-5855 (2016).

#### 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