

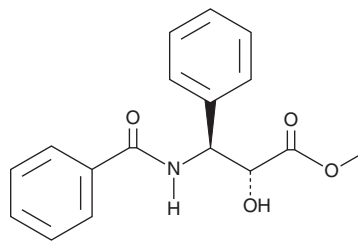
PRODUCT INFORMATION



Bz-RS-ISer(3-Ph)-OMe

Item No. 40291

CAS Registry No.: 32981-85-4
Formal Name: βS-(benzoylamino)-αR-hydroxy-benzenepropanoic acid, methyl ester
MF: C₁₇H₁₇NO₄
FW: 299.3
Purity: ≥98%
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

Bz-RS-ISer(3-Ph)-OMe is supplied as a solid. A stock solution may be made by dissolving the Bz-RS-ISer(3-Ph)-OMe in the solvent of choice, which should be purged with an inert gas. Bz-RS-ISer(3-Ph)-OMe is soluble in DMSO, acetonitrile, and methanol.

Description

Bz-RS-ISer(3-Ph)-OMe is an intermediate in the synthesis of 7-*epi*-10-deacetyltaxol (Item No. 34296), a taxane diterpenoid with α-glucosidase inhibitory and anticancer activities.¹

Reference

1. Denis, J.-N., Greene, A.E., Serra, A.A., *et al.* An efficient, enantioselective synthesis of the taxol side chain. *The Journal of Organic Chemistry* **51**(1), 46–50 (1986).

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, 03/15/2024

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