

PRODUCT INFORMATION



TP-808

Item No. 24181

CAS Registry No.: 852821-06-8
Formal Name: (4aS,8aS,9S)-9-(dimethylamino)-4a-[[[(1,1-dimethylethyl)dimethylsilyl]oxy]-8a,9-dihydro-3-(phenylmethoxy)-naphth[2,3-d]isoxazole-4,5(4aH,8H)-dione

MF: C₂₆H₃₄N₂O₅Si

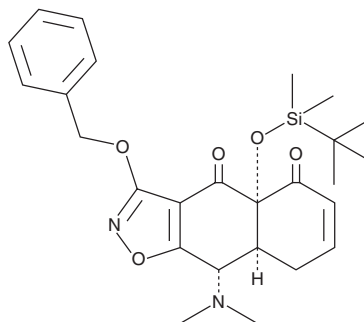
FW: 482.6

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

TP-808 is supplied as a crystalline solid. A stock solution may be made by dissolving the TP-808 in the solvent of choice, which should be purged with an inert gas. TP-808 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 20 mg/ml.

Description

TP-808 is a key intermediate in the synthesis of synthetic tetracycline antibiotics.¹

Reference

1. Zhang, W.-Y., Chen, C.-L., He, M., *et al.* Process research and development of TP-808: A key intermediate for the manufacture of synthetic tetracyclines. *Org. Process Res. Dev.* **21**(3), 377-386 (2017).

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, 11/16/2022

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