

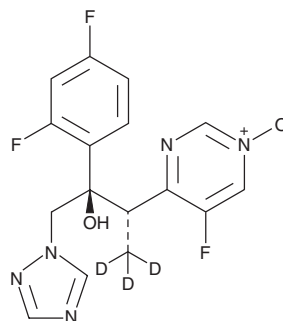
PRODUCT INFORMATION



Voriconazole-d₃ N-oxide

Item No. 40332

CAS Registry No.: 1217851-84-7
Formal Name: (αR,βS)-α-(2,4-difluorophenyl)-5-fluoro-β-methyl-α-(1H-1,2,4-triazol-1-ylmethyl)-4-pyrimidineethanol-d₃, 1-oxide
MF: C₁₆H₁₁D₃F₃N₅O₂
FW: 368.3
Chemical Purity: ≥95% (Voriconazole N-oxide)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A 1 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Voriconazole-d₃ N-oxide is intended for use as an internal standard for the quantification of voriconazole N-oxide (Item No. 28047) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Voriconazole-d₃ N-oxide is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methanol purged with an inert gas can be used. Voriconazole-d₃ N-oxide is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Voriconazole N-oxide is a major metabolite of the triazole antifungal voriconazole (Item No. 15633).¹ It is formed via oxidation of voriconazole by the cytochrome P450 (CYP) isoforms CYP3A4 and CYP2C19.

Reference

1. Hyland, R., Jones, B.C., and Smith, D.A. Identification of the cytochrome P450 enzymes involved in the N-oxidation of voriconazole. *Drug Metab. Dispos.* **31**(5), 540-547 (2003).

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, 03/25/2026

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