

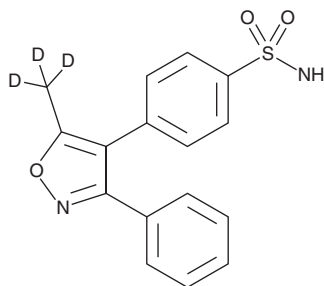
PRODUCT INFORMATION



Valdecoxib-d₃

Item No. 28699

CAS Registry No.: 1219794-90-7
Formal Name: 4-(5-(methyl-d₃)-3-phenylisoxazol-4-yl) benzenesulfonamide
MF: C₁₆H₁₁D₃N₂O₃S
FW: 317.4
Chemical Purity: ≥98% (Valdecoxib)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
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

Valdecoxib-d₃ is intended for use as an internal standard for the quantification of valdecoxib (Item No. 10006120) 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).

Valdecoxib-d₃ is supplied as a solid. A stock solution may be made by dissolving the valdecoxib-d₃ in the solvent of choice, which should be purged with an inert gas. Valdecoxib-d₃ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of valdecoxib-d₃ in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

Valdecoxib is a selective inhibitor of COX-2 (IC₅₀s = 0.005 and 140 μM, for human recombinant COX-2 and COX-1, respectively).¹ It decreases LPS-induced production of prostaglandin E₂ (PGE₂; Item No. 14010) in isolated human whole blood (IC₅₀ = 0.89 μM). Valdecoxib reduces carrageenan-induced paw edema and adjuvant-induced arthritis in rats (ED₅₀s = 10.2 and 0.032 mg/kg, respectively).

Reference

1. Talley, J.J., Brown, D.L., Carter, J.S., *et al.* 4-[5-Methyl-3-phenylisoxazol-4-yl]-benzenesulfonamide, valdecoxib: A potent and selective inhibitor of COX-2. *J. Med. Chem.* **43**(5), 775-777 (2000).

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, 01/23/2026

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