

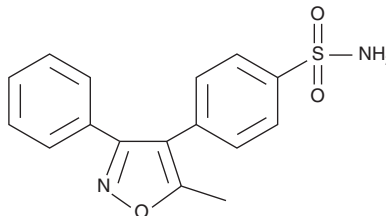
PRODUCT INFORMATION



Valdecoxib

Item No. 10006120

CAS Registry No.: 181695-72-7
Formal Name: 4-(5-methyl-3-phenyl-4-isoxazolyl)-benzenesulfonamide
Synonym: SC-65872
MF: C₁₆H₁₄N₂O₃S
FW: 314.4
Purity: ≥98%
UV/Vis.: λ_{max}: 202, 235 nm
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

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

Valdecoxib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, valdecoxib should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Valdecoxib has a solubility of approximately 0.5 mg/ml in a 1:8 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

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, 11/09/2023

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