

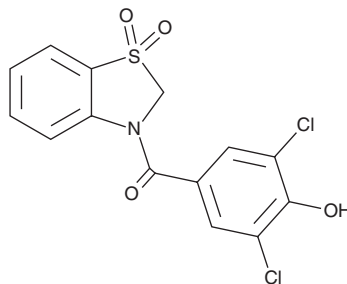
PRODUCT INFORMATION



Dotinurad

Item No. 41014

CAS Registry No.: 1285572-51-1
Formal Name: (3,5-dichloro-4-hydroxyphenyl)(1,1-dioxido-3(2H)-benzothiazolyl)-methanone
Synonym: FYU-981
MF: C₁₄H₉Cl₂NO₄S
FW: 358.2
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

Dotinurad is supplied as a solid. A stock solution may be made by dissolving the dotinurad in the solvent of choice, which should be purged with an inert gas. Dotinurad is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Dotinurad is a urate transporter 1 (URAT1) antagonist (IC₅₀ = 0.0372 μM).¹ It is selective for URAT1 over ATP-binding cassette subfamily G member 2 (ABCG2), organic anion transporter 1 (OAT1), and OAT3 (IC₅₀s = 4.16, 4.08, and 1.32 μM, respectively). *In vivo*, dotinurad (30 mg/kg) decreases plasma levels of urate and increases urine levels of urate in *Cebus* monkeys. Formulations containing dotinurad have been used in the treatment of hyperuricemia and gout.

Reference

1. Taniguchi, T., Ashizawa, N., Matsumoto, K., *et al.* Pharmacological evaluation of dotinurad, a selective urate reabsorption inhibitor. *J. Pharmacol. Exp. Ther.* **371**(1), 162-170 (2019).

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, 07/03/2024

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