

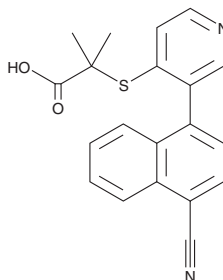
PRODUCT INFORMATION



Verinurad

Item No. 29667

CAS Registry No.: 1352792-74-5
Formal Name: 2-[[3-(4-cyano-1-naphthalenyl)-4-pyridinyl]thio]-2-methylpropanoic acid
MF: C₂₀H₁₆N₂O₂S
FW: 348.4
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

Verinurad is supplied as a crystalline solid. A stock solution may be made by dissolving the verinurad in the solvent of choice, which should be purged with an inert gas. Verinurad is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of verinurad in these solvents is approximately 30 mg/ml.

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

Description

Verinurad is an inhibitor of urate anion transporter 1 (URAT1; IC₅₀ = 0.025 μM for the human transporter).¹ It is selective for human URAT1 over rat URAT1 (IC₅₀ = 41 μM), as well as human organic anion transporter 1 (OAT1) and OAT4 (IC₅₀s = 4.6 and 5.9 μM, respectively).

Reference

1. Tan, P.K., Liu, S., Gunic, E., *et al.* Discovery and characterization of verinurad, a potent and specific inhibitor of URAT1 for the treatment of hyperuricemia and gout. *Sci. Rep.* **7**(1), 665 (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, 10/04/2022

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