

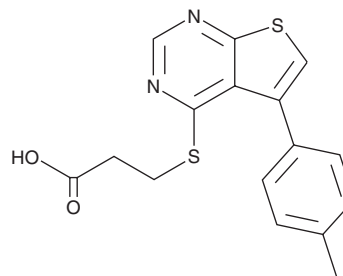
PRODUCT INFORMATION



TTP 22

Item No. 17702

CAS Registry No.: 329907-28-0
Formal Name: 3-[[5-(4-methylphenyl)thieno[2,3-d]pyrimidin-4-yl]thio]-propanoic acid
MF: C₁₆H₁₄N₂O₂S₂
FW: 330.4
Purity: ≥95%
UV/Vis.: λ_{max}: 232, 300 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

TTP 22 is supplied as a crystalline solid. A stock solution may be made by dissolving the TTP 22 in the solvent of choice, which should be purged with an inert gas. TTP 22 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TTP 22 in ethanol is approximately 2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

TTP 22 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TTP 22 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. TTP 22 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

TTP 22 is a thienopyrimidine compound that inhibits casein kinase 2 (CK2) with an IC₅₀ value of 100 nM (K_i = 40 nM).¹ It displays selectivity for CK2 over JNK3, ROCK1, and MET with no inhibitory effects towards these kinases at 10 μM.¹

Reference

1. Golub, A.G., Bdezhola, V.G., Briukhovestka, N.V., *et al.* Synthesis and biological evaluation of substituted (thieno[2,3-d]pyrimidin-4-ylthio)carboxylic acids as inhibitors of human protein kinase CK2. *Eur. J. Med. Chem.* **46**(3), 870-876 (2011).

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, 12/15/2022

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