

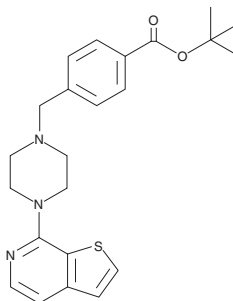
PRODUCT INFORMATION



CAY10494

Item No. 10007126

CAS Registry No.: 1044764-15-9
Formal Name: 4-[[4-thieno[2,3-c]pyridin-7-yl-1-piperazinyl)methyl]-benzoic acid, 1,1-dimethylethyl ester
MF: C₂₃H₂₇N₃O₂S
FW: 409.6
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 316 nm
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CAY10494 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10494 in the solvent of choice, which should be purged with an inert gas. CAY10494 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10494 in these solvents is approximately 5, 10, and 15 mg/ml, respectively.

CAY10494 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10494 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10494 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10494 is a synthetic intermediate useful for pharmaceutical synthesis.

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, 12/11/2022

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