

PRODUCT INFORMATION

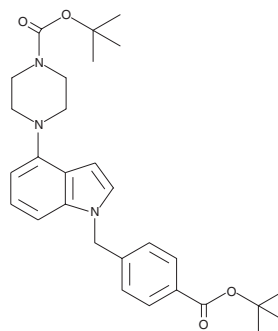


CAY10493

Item No. 10007078

CAS Registry No.: 1044764-11-5
Formal Name: 4-[1-[[4-[(1,1-dimethylethoxy) carbonyl]phenyl]methyl]-1H-indol-4-yl]-1-piperazinecarboxylic acid, 1,1-dimethylethyl ester

MF: C₂₉H₃₇N₃O₄
FW: 491.6
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 278 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

CAY10493 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10493 in the solvent of choice. CAY10493 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CAY10493 in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

CAY10493 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10493 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10493 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

CAY10493 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, 12/10/2022

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