

PRODUCT INFORMATION

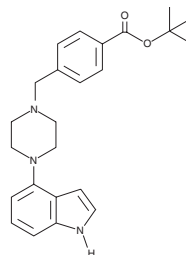


CAY10492

Item No. 10007077

CAS Registry No.: 252978-89-5
Formal Name: 4-(1H-indol-4-yl)-1-piperazine-1-carboxylic acid *tert*-butyl ester

MF: C₁₇H₂₃N₃O₂
FW: 301.4
Purity: ≥95%
UV/Vis.: λ_{max}: 222, 277 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

CAY10492 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10492 in the solvent of choice, which should be purged with an inert gas. CAY10492 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10492 in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

CAY10492 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10492 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10492 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

CAY10492 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