

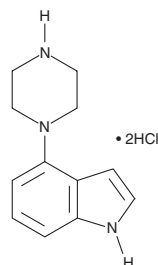
PRODUCT INFORMATION



4-(1-piperazinyl)-1H-Indole (hydrochloride)

Item No. 10007075

CAS Registry No.: 255714-24-0
Formal Name: 4-(1-piperazinyl)-1H-indole, dihydrochloride
MF: C₁₂H₁₅N₃ • 2HCl
FW: 274.2
Purity: ≥98%
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

4-(1-piperazinyl)-1H-Indole (hydrochloride) is supplied as a crystalline solid. 4-(1-piperazinyl)-1H-Indole (hydrochloride) is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of 4-(1-piperazinyl)-1H-Indole (hydrochloride) be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of 4-(1-piperazinyl)-1H-Indole (hydrochloride) in PBS, pH 7.2, is at least 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4-(1-piperazinyl)-1H-Indole (hydrochloride) 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