

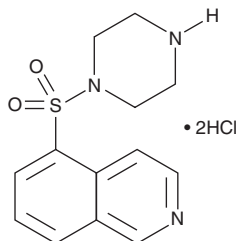
PRODUCT INFORMATION



HA-100 (hydrochloride)

Item No. 14955

CAS Registry No.: 210297-47-5
Formal Name: 5-(1-piperazinylsulfonyl)-isoquinoline, dihydrochloride
Synonym: C-1
MF: C₁₃H₁₅N₃O₂S • 2HCl
FW: 350.3
Purity: ≥95%
UV/Vis.: λ_{max}: 216, 275, 323 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

HA-100 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the HA-100 (hydrochloride) in the solvent of choice. HA-100 (hydrochloride) is soluble in DMSO at a concentration of approximately 1.1 mg/ml.

HA-100 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HA-100 (hydrochloride) should first be dissolved in dimethyl formamide (DMF) and then diluted with the aqueous buffer of choice. HA-100 (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HA-100 is an isoquinoline compound with an added piperazinylsulfonyl group that acts as an inhibitor of protein kinases (PKs), including PKA, PKC, and PKG (IC₅₀s = 8, 12, and 4 μM, respectively).¹ It less effectively blocks the activity of myosin light chain kinase (IC₅₀ = 240 μM).¹

Reference

1. Hagiwara, M., Inagaki, M., Watanabe, M., *et al.* Selective modulation of calcium-dependent myosin phosphorylation by novel protein kinase inhibitors, isoquinolinesulfonamide derivatives. *Mol. Pharmacol.* **32**(1), 7-12 (1987).

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, 11/15/2022

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