

PRODUCT INFORMATION



KN-92 (hydrochloride)

Item No. 9000890

CAS Registry No.: 1431698-47-3
Formal Name: 2-[N-(4-methoxybenzenesulfonyl)]

amino-N-(4-chlorocinnamyl)-
N-methylbenzylamine,
monohydrochloride

MF: $C_{24}H_{25}ClN_2O_3S \cdot HCl$

FW: 493.4

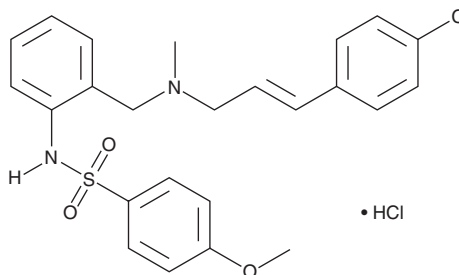
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 204, 251 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

KN-92 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the KN-92 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. KN-92 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KN-92 (hydrochloride) in these solvents is approximately 10 mg/ml.

Description

KN-92 is an inactive derivative of KN-93, the selective inhibitor of Ca^{2+} /calmodulin-dependent kinase type II (CaMKII).¹ At concentrations up to 25 μM , KN-92 is ineffective at inhibiting CaMKII or arresting cell growth of NIH 3T3 fibroblasts.² It is intended to be used as a control compound in studies designed to elucidate the antagonist activities of KN-93 (Item No. 13319).

References

- Sumi, M., Kiuchi, K., Ishikawa, T., *et al.* The newly synthesized selective Ca^{2+} /calmodulin dependent protein kinase II inhibitor KN-93 reduces dopamine contents in PC12h cells. *Biophys. Res. Commun.* **181**(3), 968-975 (1991).
- Tombes, R.M., Grant, S., Westin, E.H., *et al.* G1 cell cycle arrest and apoptosis are induced in NIH 3T3 cells by KN-93, an inhibitor of CaMK-II (the multifunctional Ca^{2+} /CaM kinase). *Cell Growth Differ.* **6**(9), 1063-70 (1995).

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, 04/26/2024

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