

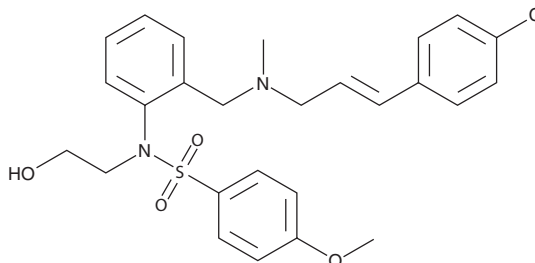
Product Information



KN-93

Catalog No. 13319

CAS Registry No.: 139298-40-1
Formal Name: N-[2-[[[3-(4-chlorophenyl)-2-propen-1-yl]methylamino]methyl]phenyl]-N-(2-hydroxyethyl)-4-methoxybenzenesulfonamide
MF: C₂₆H₂₉ClN₂O₄S
FW: 501.0
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that KN-93 be stored as supplied at -20°C. It should be stable for at least two years.

KN-93 is supplied as a crystalline solid. A stock solution may be made by dissolving the KN-93 in an organic solvent purged with an inert gas. KN-93 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of KN-93 in these solvents is approximately 30 mg/ml.

KN-93 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, KN-93 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. KN-93 has a solubility of approximately 0.25 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

KN-93 is a water-soluble, selective inhibitor of Ca²⁺/calmodulin-dependent kinase II (CaMKII), competitively blocking CaM binding to the kinase (K_i = 370 nM).¹ It does not affect the activities of PKA, PKC, MLCK, or Ca²⁺-phosphodiesterase.¹ It inhibits histamine-induced aminopyrine uptake in parietal cells (IC₅₀ = 300 nM).² More recently, KN-93 has been used to implicate roles for CaMKII in Ca²⁺-induced Ca²⁺ release in cardiac myocytes, constitutive phosphorylation of 5-lipoxygenase in 3T3 cells, and Ca²⁺-dependent activation of HIF-1α in colon cancer cells.³⁻⁵

References

- Sumi, M., Kiuchi, K., Ishikawa, T., *et al.* The newly synthesized selective Ca²⁺/calmodulin dependent protein kinase II inhibitor KN-93 reduces dopamine contents in PC12h cells. *Biochemical and Biophysical Research Communications* **181**(3), 968-975 (1991).
- Mamiya, N., Goldenring, J.R., Tsunoda, Y., *et al.* Inhibition of acid secretion in gastric parietal cells by the Ca²⁺/calmodulin-dependent protein kinase II inhibitor KN-93. *Biochemical and Biophysical Research Communications* **195**(2), 608-615 (1993).
- Oestreich, E.A., Malik, S., Goonasekera, S.A., *et al.* Epac and phospholipase Cε regulate Ca²⁺ release in the heart by activation of protein kinase Cε and calcium-calmodulin kinase II. *J. Biol. Chem.* **284**(3), 1514-1522 (2009).
- Flamand, N., Luo, M., Peters-Golden, M., *et al.* Phosphorylation of serine 271 on 5-lipoxygenase and its role in nuclear export. *J. Biol. Chem.* **284**(1), 306-313 (2009).
- Riganti, C., Doublier, S., Viarisio, D., *et al.* Artemisinin induces doxorubicin resistance in human colon cancer cells via calcium-dependent activation of HIF-1α and P-glycoprotein overexpression. *Brit. J. Pharmacol.* **156**, 1054-1066 (2009).

Related Product

KN-62 - Cat. No. 13318

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any **obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 09/03/2009

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com