

# PRODUCT INFORMATION



## KN-93 (phosphate)

Item No. 21472

**CAS Registry No.:** 1188890-41-6  
**Formal Name:** N-[2-[[[(2E)-3-(4-chlorophenyl)-2-propen-1-yl]methylamino]methyl]phenyl]-N-(2-hydroxyethyl)-4-methoxy-benzenesulfonamide, monophosphate

**MF:** C<sub>26</sub>H<sub>29</sub>ClN<sub>2</sub>O<sub>4</sub>S • H<sub>3</sub>PO<sub>4</sub>

**FW:** 599.0

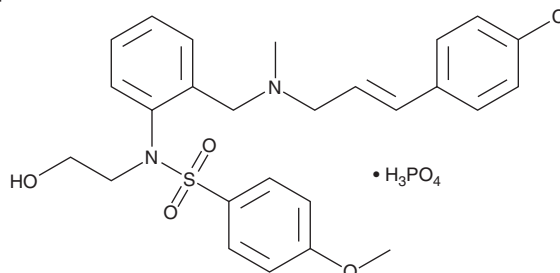
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 252 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

KN-93 (phosphate) is supplied as a crystalline solid. A stock solution may be made by dissolving the KN-93 (phosphate) in the solvent of choice, which should be purged with an inert gas. KN-93 (phosphate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KN-93 (phosphate) in these solvents is approximately 30 and 25 mg/ml, respectively. KN-93 (phosphate) is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of KN-93 (phosphate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. KN-93 (phosphate) is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

KN-93 (phosphate) is a potent and selective inhibitor of Ca<sup>2+</sup>/calmodulin-dependent kinase II (CaMKII) (K<sub>i</sub> = 370 nM), inhibiting both the α- and β-subunits of CaMKII.<sup>1</sup> It does not have significant effects on cAMP-dependent protein kinase, Ca<sup>2+</sup>/phospholipid-dependent protein kinase, myosin light chain kinase, or Ca<sup>2+</sup> phosphodiesterase activity. It inhibits histamine-induced aminopyrine uptake in parietal cells (IC<sub>50</sub> = 300 nM).<sup>2</sup> KN-93 can block voltage-gated potassium (K<sub>v</sub>) channels when applied extracellularly, independent of its CaMKII action.<sup>3</sup> More recently, KN-93 has been used to implicate roles for CaMKII in Ca<sup>2+</sup>-induced Ca<sup>2+</sup> release in cardiac myocytes, constitutive phosphorylation of 5-lipoxygenase in 3T3 cells, and Ca<sup>2+</sup>-dependent activation of HIF-1α in colon cancer cells.<sup>4-6</sup>

### References

1. Sumi, M., Kiuchi, K., Ishikawa, T., *et al.* *Biochem. Biophys. Res. Commun.* **181(3)**, 968-975 (1991).
2. Mamiya, N., Goldenring, J.R., Tsunoda, Y., *et al.* *Biochem. Biophys. Res. Commun.* **195(2)**, 608-615 (1993).
3. Rezazadeh, S., Claydon, T.W. and Fedida, D.J. *Pharmacol. Exp. Ther.* **317(1)**, 292-299 (2006).
4. Oestreich, E.A., Malik, S., Goonasekera, S.A., *et al.* *J. Biol. Chem.* **284(3)**, 1514-1522 (2009).
5. Flamand, N., Luo, M., Peters-Golden, M., *et al.* *J. Biol. Chem.* **284(1)**, 306-313 (2009).
6. Riganti, C., Doublie, S., Viariso, D., *et al.* *Br. J. Pharmacol.* **156(7)**, 1054-1066 (2009).

#### 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, 11/23/2022

#### 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