# **KN-93**

#### **CaM** kinase inhibitor

KN-93 is a potent, selective and cell-permeable inhibitor of CaM kinase II (IC $_{50}$ =370 nM). Does not significantly inhibit PKA. Causes cell-cycle arrest in NIH 3T3 fibroblasts and HeLa cells. Inhibits glucose-induced and forskolin-stimulated insulin release from isolated pancreatic islets of Langerhans. Induces  $\rm G_1$  cell cycle arrest and apoptosis in NIH 3T3 cells.

Citations: 9

View Online »

**Ordering Information** 

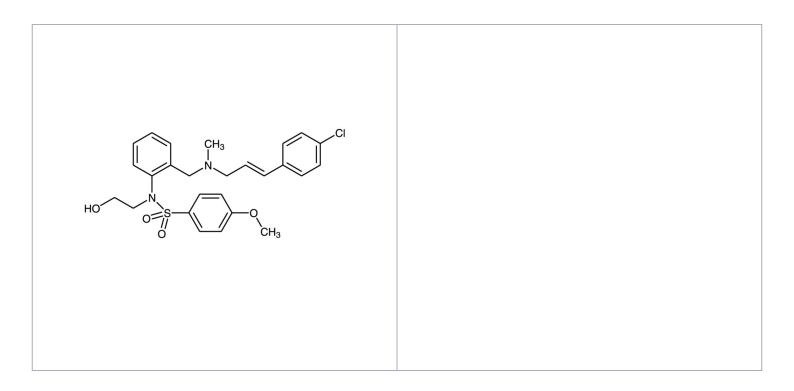
Order Online »

BML-EI268-0001

1mg

Manuals, SDS & CofA

**View Online** »



### **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended. Stable for at least

2 years after receipt. Stock solutions are stale for up to 3 months when stored at -20°C.

Handling Protect from light. Hygroscopic.

**Long Term Storage** -20°C

**Shipping Ambient Temperature** 

## Regulatory Status RUO - Research Use Only

#### **Product Details**

**Appearance** white to off-white solid

**CAS** 139298-40-1

**Couple Target** CaM kinase

**Couple Type** Inhibitor

 $C_{26}H_{29}CIN_2O_4S$ **Formula** 

MW 501

**Purity** ≥95% (HPLC)

**Solubility** Soluble in DMSO (25mg/ml).

Last modified: May 29, 2024

European Sales Office