# K-252a

#### Protein kinase inhibitor

Alkaloid isolated from soil fungi. General, cell permeable protein kinase inhibitor. Acts by binding to the ATP binding domain of the kinase. Potent inhibitor of  ${\rm Ca^{2^+}/calmodulin}$  kinase II. Inhibits myosin light chain kinase, cAMP-dependent protein kinase (PKA), protein kinase C (PKC) (K $_{\rm i}$ =25nM), trk tyrosine kinase family members (gp140trk; IC $_{\rm 50}$  =3nM) and cGMP-dependent protein kinase (PKG). Binds with high affinity to neuroblastoma membrane (2.7nM), promoting survival of primary neuronal cultures (75nM) and induction of neurite outgrowth in SH-SY5Y neuroblastoma cells along with phosphorylation of pp125 focal adhesion protein tyrosine kinase. Induces apoptosis.

Citations: 30

View Online »

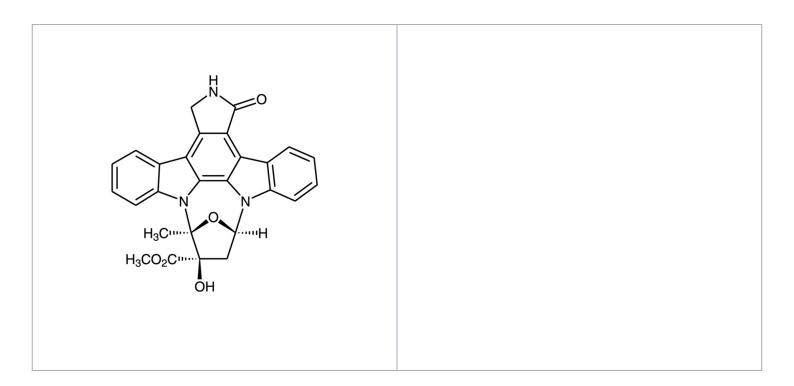
## **Ordering Information**

Order Online »

BML-EI152-1000	1mg
BML-EI152-0100	100µg

Manuals, SDS & CofA

**View Online »** 



## **Handling & Storage**

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

1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up

to 3 months at -20°C.

**Handling** Protect from light.

Long Term Storage -20°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

#### **Product Details**

**Appearance** Lyophilized solid.

**CAS** 99533-80-9

Couple Target CaM kinase, MLCK, PKA, PKC, PKG, Trk

Couple Type Inhibitor

Formula  $C_{27}H_{21}N_3O_5$ 

**MW** 467.5

Purity ≥98% (HPLC)

RTECS NZ0550000

Soluble in DMSO (1mg/ml) or dimethyl

formamide (1mg/ml).

Last modified: May 29, 2024

