

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 Ca^{2+} /calmodulin kinase II. Inhibits myosin light chain kinase, cAMP-dependent protein kinase (PKA), protein kinase C (PKC) ($K_i=25\text{nM}$), trk tyrosine kinase family members (gp140trk; $\text{IC}_{50}=3\text{nM}$) 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 »](#)

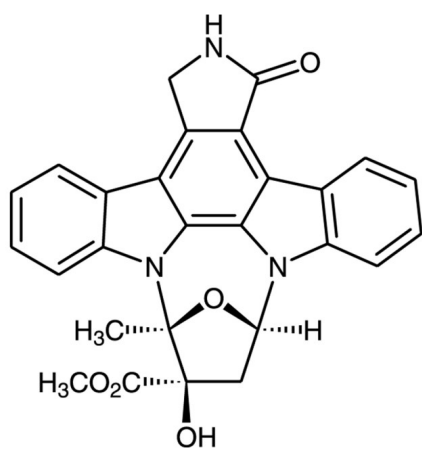
Ordering Information

[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
Solubility	Soluble in DMSO (1mg/ml) or dimethyl formamide (1mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)