

GF109203X

PKC inhibitor

Potent and selective protein kinase C inhibitor which has been used effectively in platelets, Swiss 3T3 fibroblasts and macrophages. Inhibitory profile: PKC, $IC_{50}=0.02\mu M$ (inhibits α , βI , βII and γ subtypes with similar potency); PKA, $IC_{50}=2.0\mu M$; phosphorylase kinase, $IC_{50}=0.7\mu M$; insulin, EGF and PDGF receptor tyrosine kinases are not inhibited at concentrations up to $50\mu M$.

Citations: 21

[View Online »](#)

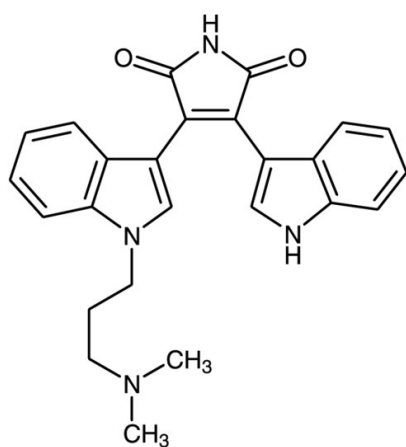
Ordering Information

[Order Online »](#)

BML-EI246-0001	1mg
BML-EI246-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-[1-(3-Dimethylaminopropyl)-1H-indol-3-yl]-3-(1H-indol-3-yl)maleimide, Bisindolylmaleimide I
Appearance	Red solid.
CAS	133052-90-1
Couple Target	PKC
Couple Type	Inhibitor
Formula	$C_{25}H_{24}N_4O_2$
MW	412.5
Purity	≥98% (HPLC)
Solubility	Soluble in dimethyl formamide or DMSO (10mg/ml, warm to 40°C and vortex well).

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)