

KT5823

PKG inhibitor

KT5823 is a derivative of K-252a and is a selective, cell permeable, inhibitor of PKG ($K_i=0.234 \mu\text{M}$). K_i 's for PKA = $4\mu\text{M}$ and for MLCK => $10\mu\text{M}$.

Citations: 3

[View Online »](#)

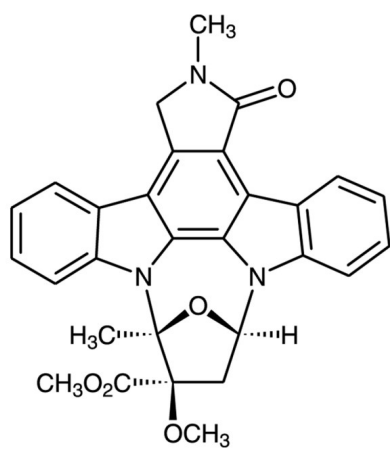
Ordering Information

[Order Online »](#)

BML-EI200-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at 4°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	126643-37-6
Couple Target	PKG
Couple Type	Inhibitor
Formula	$C_{29}H_{25}N_3O_5$
MW	495.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (20mg/ml), dimethyl formamide (20mg/ml) and ethyl acetate (5mg/ml).

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)