

# 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 »](#)

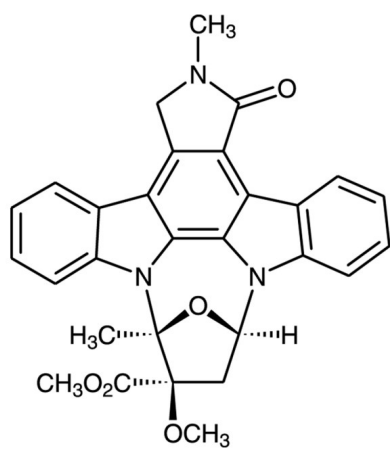
## Ordering Information

[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).



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)