

# Go-6983

## PKC inhibitor

PKC inhibitor. Has been shown to selectively inhibit several PKC isoenzymes ( $IC_{50}$  = 7nM for PKC $\alpha$  and PKC $\beta$ ; 6nM for PKC $\gamma$ ; 10nM for PKC $\delta$ ; 60nM for PKC $\zeta$ ). The compound does not effectively inhibit PKC $\mu$  ( $IC_{50}$ =20  $\mu$ M) and therefore can be used to differentiate PKC $\mu$  from other isoforms.

Citations: 3

[View Online »](#)

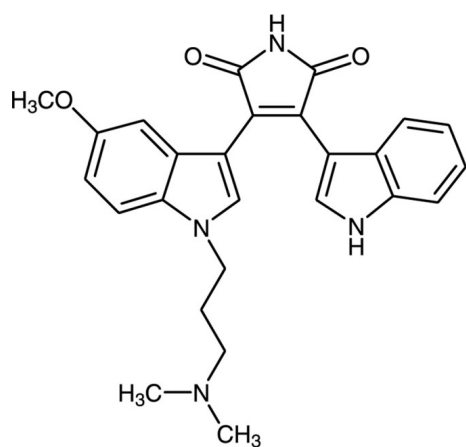
## Ordering Information

[Order Online »](#)

BML-EI296-0500	0.5mg
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** Bright red solid.

**CAS** 133053-19-7

**Couple Target** PKC

**Couple Type** Inhibitor

**Formula**  $C_{26}H_{26}N_4O_3$

**Identity** Determined by  $^1H$ -NMR and MS.

**MW** 442.5

**Purity**  $\geq 95\%$  (HPLC)

**Solubility** Soluble in DMSO.

**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-  
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)