

# Ro 31-8220

## PKC inhibitor

Cell-permeable PKC inhibitor. Selective for PKC ( $IC_{50}=10nM$ ) over CaM kinase II ( $IC_{50}=17\mu M$ ) and PKA ( $IC_{50}=0.9\mu M$ ). Induces apoptosis in various cell lines.

Citations: 10

[View Online »](#)

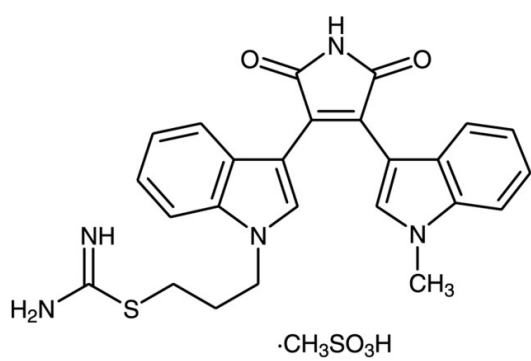
## Ordering Information

[Order Online »](#)

|                |     |
|----------------|-----|
| BML-EI283-0001 | 1mg |
|----------------|-----|

## 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-(Amidinothio)propyl)-1H-indol-3-yl]-3-(1-methylindol-3-yl)maleimide·methanesulfonate, Bisindolylmaleimide IX . Methanesulfonate |
| Appearance       | Red/orange solid.                                                                                                                       |
| CAS              | 138489-18-6                                                                                                                             |
| Couple Target    | PKC                                                                                                                                     |
| Couple Type      | Inhibitor                                                                                                                               |
| Formula          | $C_{25}H_{23}N_5O_2S \cdot CH_4O_3S$                                                                                                    |
| MW               | 553.7                                                                                                                                   |
| Purity           | ≥98%                                                                                                                                    |
| Solubility       | Soluble in DMSO (10mg/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-be@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)