

# Chelerythrine chloride

PKC inhibitor

Reported to be a potent and selective inhibitor of protein kinase C ( $IC_{50} = 0.66\mu M$ ). Although chelerythrine has numerous documented biological effects, studies have raised serious doubts about the effectiveness of chelerythrine as a PKC inhibitor.

Citations: 10

[View Online »](#)

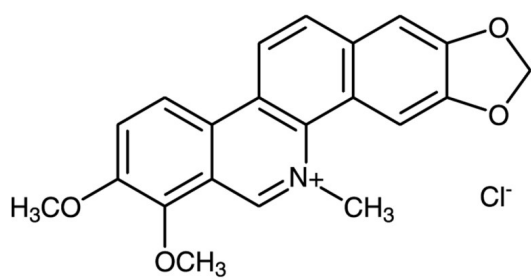
## Ordering Information

[Order Online »](#)

BML-EI225-0005	5mg
BML-EI225-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Heating to 40-50°C with gentle agitation may support complete solubilization. Further dilute with buffer just prior to use. Stock solutions should be refrigerated or frozen between for storage.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Yellow solid.
CAS	3895-92-9
Couple Target	PKC
Couple Type	Inhibitor
Formula	$C_{21}H_{18}ClNO_4$
MI	14: 2051
MW	383.8
Purity	≥98% (HPLC, TLC)
RTECS	FL9200000
Solubility	Soluble in DMSO, water (2mg/ml warm), 100% ethanol or methanol.

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-  
be@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)