

# Bryostatin 1

## PKC activator

A structurally unique marine natural product that binds to and activates PKC at picomolar concentrations, but is not a carcinogen or a complete tumor promoter. In many systems it induces only a subset of the responses to PMA and blocks those which it does not induce. It is markedly more potent than PMA for translocating PKC $\delta$  and  $\epsilon$ .

Citations: 6

[View Online »](#)

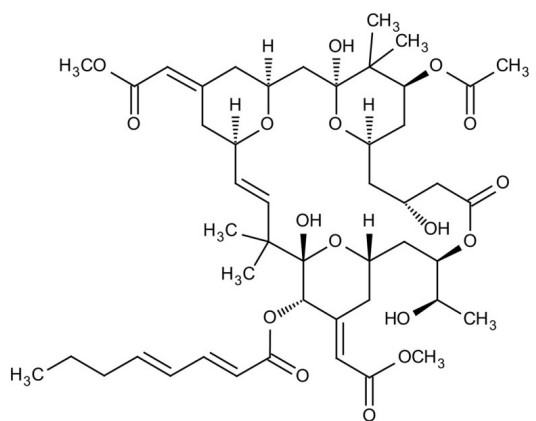
## Ordering Information

[Order Online »](#)

BML-ST103-0010	10 $\mu$ g
----------------	------------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	White solid to colorless film.
CAS	83314-01-6
Couple Target	PKC
Couple Type	Activator
Formula	$C_{47}H_{68}O_{17}$
MW	905.0
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/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-  
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)