

# D609

## Phospholipase inhibitor

Selective inhibitor of phosphatidylcholine-specific phospholipase C (PC-PLC). Has been used to study the coupling of phosphatidylcholine-specific phospholipase C with sphingomyelinase. Shows antitumor and antiviral activity. Inhibits basement membrane collagen synthesis, which is the last step in the development of a new blood vessel. Inhibits induction of nitric oxide synthases (NOS). Induces apoptosis. It indirectly inhibits phorbol ester stimulated PLD activity.

Citations: 26

[View Online »](#)

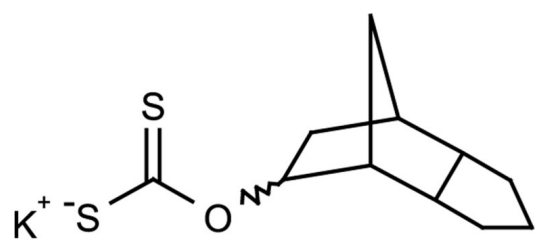
## Ordering Information

[Order Online »](#)

BML-ST330-0005	5mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. We recommend to prepare fresh solutions each day.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Tricyclodecan-9-yl xanthogenate . K
<b>Appearance</b>	Off-white solid.
<b>CAS</b>	83373-60-8
<b>Couple Target</b>	Phospholipase, PLC
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{11}H_{15}OS_2 \cdot K$
<b>MW</b>	266.5
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in water (25mg/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-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)