

# SQ 22536

## Adenylate cyclase inhibitor

SQ 22536 is a cell-permeable adenylate cyclase inhibitor. It inhibits cAMP production induced by prostaglandin E<sub>1</sub> in platelet lysates (IC<sub>50</sub>=82 μM) and in intact platelets (IC<sub>50</sub>=1 μM). It has been successfully employed as a probe for the involvement of adenylate cyclase in a variety of receptor systems including those for prostaglandins and neurotransmitters. Reduces PGE<sub>2</sub>-induced inhibition of O<sub>2</sub><sup>-</sup> production.

Citations: 18

[View Online »](#)

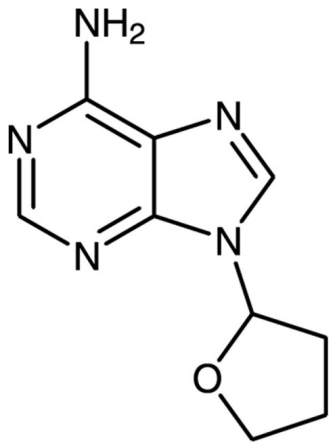
## Ordering Information

[Order Online »](#)

BML-CN140-0005	5mg
BML-CN140-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	9-THF-Ade
<b>Appearance</b>	White to off-white solid.
<b>CAS</b>	17318-31-9
<b>Couple Target</b>	Adenylate cyclase
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_9H_{11}N_5O$
<b>MW</b>	205.2
<b>Purity</b>	≥99% (TLC)
<b>Solubility</b>	Soluble in water (25mg/ml) or DMSO (50mM with warming).
<b>Source</b>	Synthetic.



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)