

SQ 22536

Adenylate cyclase inhibitor

SQ 22536 is a cell-permeable adenylate cyclase inhibitor. It inhibits cAMP production induced by prostaglandin E₁ in platelet lysates (IC₅₀=82 μM) and in intact platelets (IC₅₀=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 PGE₂-induced inhibition of O₂⁻ production.

Citations: 18

[View Online »](#)

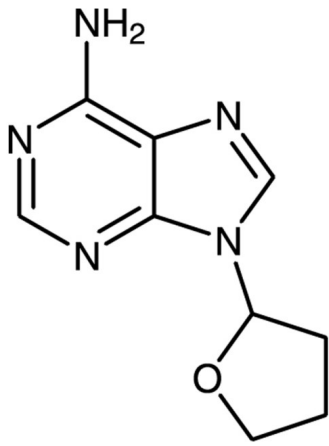
Ordering Information

[Order Online »](#)

BML-CN140-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status RUO - Research Use Only

Product Details

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com