

# Adenosine 5'-O-(3-thiotriphosphate) . tetralithium salt

**K<sup>+</sup> channel activator**

ATP analog that can substitute ATP in various kinase reactions. Substrate and inhibitor of ATP-dependent enzyme systems. The resulting thiophosphorylated proteins are resistant to dephosphorylation by protein phosphatases. Increases the activity of Ca<sup>2+</sup>-activated K channels.

Citations: 6

[View Online »](#)

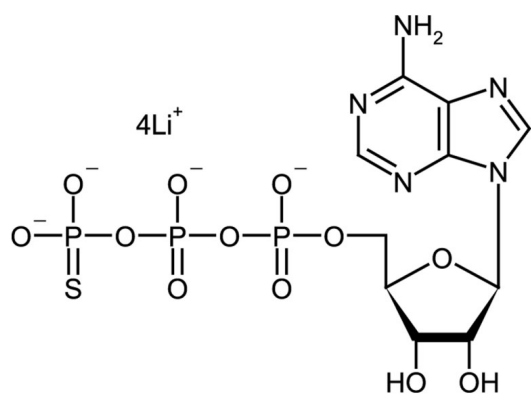
## Ordering Information

[Order Online »](#)

ALX-480-066-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions may be stored at -80°C for up to 1 month. Product is not sterile.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	ATP-γS . 4Li
Appearance	White to off-white solid.
CAS	93839-89-5
Couple Target	Potassium channel
Couple Type	Activator
Formula	$C_{10}H_{12}N_5O_{12}P_3S \cdot 4Li$
MW	547
Purity	≥90% (HPLC)
Solubility	Soluble in water.



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-  
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)