

# 2-

# Aminoethoxydiphenylborate

## TRP channel activator

Cell permeable modulator of  $\text{Ins}(1,4,5)\text{-P}_3$ -induced  $\text{Ca}^{2+}$  release. A prototype drug for a group of structurally related calcium channel blockers in human platelets. Has been shown, in the absence of other stimuli, to activate TRPV1, V2 and V3, but not TRPV4, V5 and V6 expressed in HEK 293 cells. Potentiates the effect of capsaicin, protons and heat on TRPV1. It was originally introduced as an IP3 receptor blocker and subsequently as an inhibitor of store-operated  $\text{Ca}^{2+}$  entry (75 $\mu\text{M}$ ).

Citations: 21

[View Online »](#)

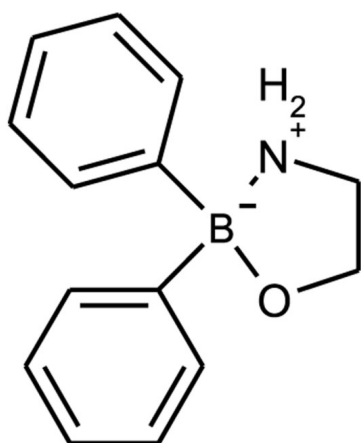
## Ordering Information

[Order Online »](#)

ALX-400-045-M100	100mg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 2 months when stored at -20°C. We do not recommend storing aqueous solutions for more than one day.
Handling	Packaged under inert gas. Hygroscopic. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2-APB
Appearance	White to off-white crystalline solid.
CAS	524-95-8
Couple Target	IP3 receptor, TRP channel
Couple Type	Activator, Inhibitor
Formula	$C_{14}H_{16}BNO$
MW	225.1
Purity	≥98% (HPLC)
RTECS	ED6150000
Solubility	Soluble in DMSO (20mg/ml), dimethyl formamide (20mg/ml) or 100% ethanol (5mg/ml); slightly soluble in PBS, pH 7.2 (50µg/ml).



ENZO LIFE SCIENCES,  
INC.  
Phone: +1 949 949 9120  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

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