

# CBA

## TRPM4 blocker

CBA is a selective blocker of TRPM4 channels with no activity at TRPM5, TRPM7, TRPM8, TRPV1, TRPV3, TRPV6 and a panel of other ion channels and receptors. Displays protective effects in murine experimental autoimmune encephalomyelitis and in a model of glutamate-induced neuronal degeneration. A highly useful tool for exploring the functions of TRPM4 in a variety of cell types.

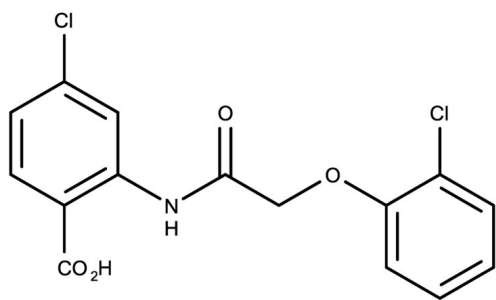
### Ordering Information

[Order Online »](#)

ENZ-CHM446-0025	25mg
-----------------	------

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 4-Chloro-2-[[2-(2-chlorophenoxy)acetyl]amino]benzoic acid

**Appearance** Off-white solid.

**CAS** 351424-20-9

**Couple Type** Inhibitor

**Formula**  $C_{15}H_{11}Cl_2NO_4$

**Identity** Determined by NMR.

**MW** 340.16

**Purity** ≥98% (TLC)

**Solubility** Soluble in DMSO (35 mg/ml).

**Last modified: January 24, 2025**



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)