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

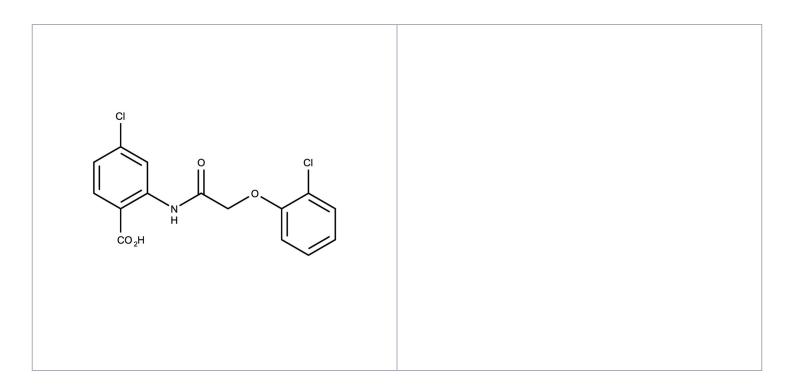
Order Online »

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 $C_{15}H_{11}CI_2NO_4$

Identity Determined by NMR.

MW 340.16

Purity ≥98% (TLC)

Solubility Soluble in DMSO (35 mg/ml).

Last modified: January 24, 2025