SK&F 96365

TRP channel blocker

A novel, selective inhibitor of receptor-mediated calcium entry. It inhibits receptor-mediated calcium entry in stimulated platelets (IC $_{50}$ =8-12µM) neutrophils and endothelial cells at concentrations which do not affect internal Ca $^{2+}$ release. Note, in some cells SK&F 96365 can activate a novel Ca $^{2+}$ entry pathway. Blocks TRPC3 and TRPC6 activation. Inhibits the histamine-induced formation of nitric oxide in human endothelial cells.

Citations: 8

View Online »

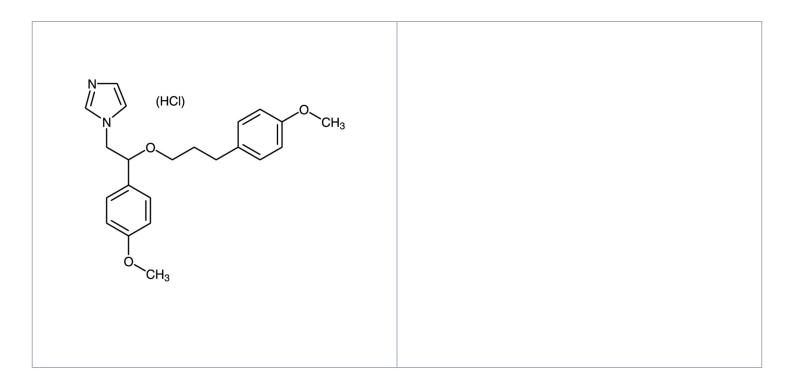
Ordering Information

Order Online »

BML-CA230-0005	5mg
BML-CA230-0025	25mg

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, as supplied, at room temperature. Stock solutions are

stable for up to 3 months at -20°C.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-[β-[3-(4-Methoxyphenyl)propoxy]-4-methoxyphenethyl]-

1H-imidazole . HCI, SKF-96365 . hydrochloride

Appearance White solid.

CAS 130495-35-1

Couple Target Calcium, TRP channel

Couple Type Blocker

Formula $C_{22}H_{26}N_2O_3$. HCI

MW 402.9

MeltingPoint 117-119°C

Purity ≥98% (HPLC)

Solubility Soluble in water (30mg/ml).

Last modified: May 29, 2024

