

# SK&F 96365

## TRP channel blocker

A novel, selective inhibitor of receptor-mediated calcium entry. It inhibits receptor-mediated calcium entry in stimulated platelets ( $IC_{50}$ =8-12 $\mu$ M) 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 »](#)

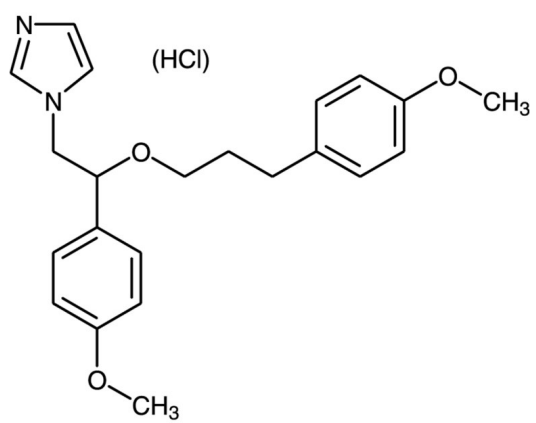
## Ordering Information

[Order Online »](#)

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

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	1-[β-[3-(4-Methoxyphenyl)propoxy]-4-methoxyphenethyl]-1H-imidazole . HCl, SKF-96365 . hydrochloride
<b>Appearance</b>	White solid.
<b>CAS</b>	130495-35-1
<b>Couple Target</b>	Calcium, TRP channel
<b>Couple Type</b>	Blocker
<b>Formula</b>	$C_{22}H_{26}N_2O_3 \cdot HCl$
<b>MW</b>	402.9
<b>MeltingPoint</b>	117-119°C
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in water (30mg/ml).



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)