

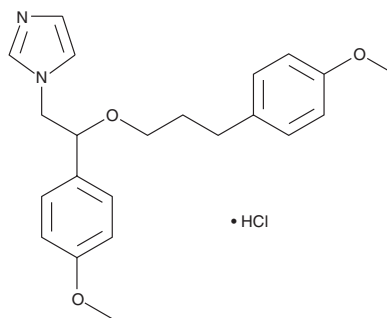
PRODUCT INFORMATION



SKF 96365 (hydrochloride)

Item No. 10009312

CAS Registry No.: 130495-35-1
Formal Name: 1-[2-(4-methoxyphenyl)-2-[3-(4-methoxyphenyl)propoxy]ethyl]-1H-imidazole, monohydrochloride
MF: C₂₂H₂₆N₂O₃ • HCl
FW: 402.9
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 275 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SKF 96365 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SKF 96365 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SKF 96365 (hydrochloride) is miscible in organic solvents such as ethanol, DMSO, and dimethyl formamide

SKF 96365 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SKF 96365 (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SKF 96365 (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SKF 96365 inhibits the receptor-mediated influx of calcium via voltage-gated calcium channels with an IC₅₀ value of approximately 10 μM.¹ It inhibits the acetylcholine-induced depolarization of circular smooth muscle in a dose-dependent manner at 3-50 μM.² SKF 96365 can distinguish receptor-mediated release in platelets and neutrophils from the calcium release from internal stores. However, it does not distinguish between receptor-mediated and voltage-gated release.

References

1. Merritt, J.E., Armstrong, W.P., Benham, C.D., *et al.* SK&F 96365, a novel inhibitor of receptor-mediated calcium entry. *Biochem J.* **271**, 515-522 (1990).
2. Hotta, A., Kim, Y.C., Kakamura, E., *et al.* Effects of inhibitors of nonselective cation channels on the acetylcholine-induced depolarization of circular smooth muscle from the guinea-pig stomach antrum. *J. Smooth Muscle Res.* **41(6)**, 313-327 (2005).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/03/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM