

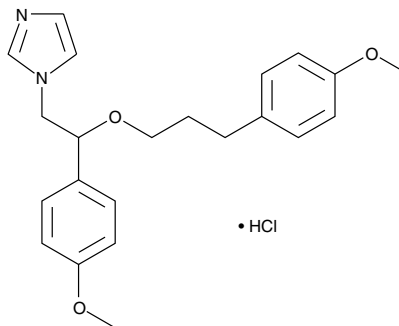
Product Information



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%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 226, 275 nm



Laboratory Procedures

For long term storage, we suggest that SKF 96365 (hydrochloride) be stored as supplied at -20°C. It should be stable for at least two years.

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

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.

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

- 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).
- 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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10009312

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 10/15/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com