

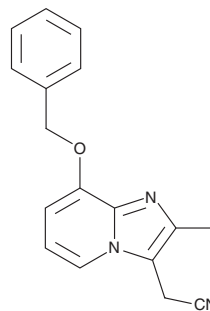
PRODUCT INFORMATION



SCH 28080

Item No. 17885

CAS Registry No.: 76081-98-6
Formal Name: 2-methyl-8-(phenylmethoxy)-imidazo[1,2-a]pyridine-3-acetonitrile
MF: C₁₇H₁₅N₃O
FW: 277.3
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 235, 280 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

SCH 28080 is supplied as a crystalline solid. A stock solution may be made by dissolving the SCH 28080 in the solvent of choice, which should be purged with an inert gas. SCH 28080 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SCH 28080 in these solvents is approximately 20 mg/ml.

SCH 28080 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SCH 28080 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SCH 28080 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

SCH 28080 is a reversible K⁺-competitive inhibitor of H⁺/K⁺-ATPases that is best known for its ability to block acid secretion through its action against the gastric H⁺/K⁺-ATPase (IC₅₀ = 1.3 μM).¹⁻³ It is effective against gastric H⁺/K⁺-ATPases from a variety of species and can inhibit colonic H⁺/K⁺-ATPases, but this activity appears to be species-dependent.⁴ SCH 28080 is also used to investigate the role of H⁺/K⁺-ATPases in non-mammalian organisms and to distinguish the actions of H⁺/K⁺-ATPases from other ATP-dependent transporters.^{5,6}

References

1. Beil, W., Hackbarth, I., and Sewing, K.-F. Mechanism of gastric antisecretory effect of SCH 28080. *Br. J. Pharmac.* **88(1)**, 19-23 (1986).
2. Wallmark, B., Briving, C., Fryklund, J., et al. Inhibition of gastric H⁺,K⁺-ATPase and acid secretion by SCH 28080, a substituted pyridyl(1,2a)imidazole. *J. Biol. Chem.* **262(5)**, 2077-2084 (1987).
3. Swarts, H.G.P., Hermesen, H.P.H., Koenderink, J.B., et al. Conformation-dependent inhibition of gastric H⁺,K⁺-ATPase by SCH 28080 demonstrated by mutagenesis of glutamic acid 820. *Mol. Pharmacol.* **55(3)**, 541-547 (1999).
4. Shao, J., Gumz, M.L., Cain, B.D., et al. Pharmacological profiles of the murine gastric and colonic H,K-ATPases. *Biochim. Biophys. Acta* **1800(9)**, 906-911 (2010).
5. Beane, W.S., Morokuma, J., Adams, D.S., et al. A chemical genetics approach reveals H,K-ATPase-mediated membrane voltage is required for planarian head regeneration. *Chem. Biol.* **18(1)**, 77-89 (2011).
6. Salyer, S.A., Olderding, J.R., Distler, A.A., et al. Vacuolar ATPase driven potassium transport in highly metastatic breast cancer cells. *Biochim. Biophys. Acta* **1832(10)**, 1734-1743 (2013).

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, 11/09/2022

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