

PRODUCT INFORMATION

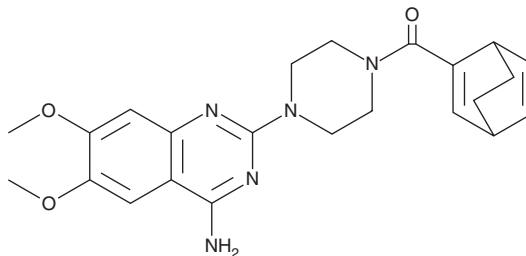


Prazobind

Item No. 21409

CAS Registry No.: 107021-36-3
Formal Name: [4-(4-amino-6,7-dimethoxy-2-quinazoliny)-1-piperazinyl] bicyclo[2.2.2]octa-2,5-dien-2-yl-methanone

Synonym: SZL-49
MF: C₂₃H₂₇N₅O₃
FW: 421.5
Purity: ≥95%
UV/Vis.: λ_{max}: 212, 247, 331 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

Prazobind is supplied as a crystalline solid. A stock solution may be made by dissolving the prazobind in the solvent of choice, which should be purged with an inert gas. Prazobind is soluble in the organic solvent DMSO at a concentration of approximately 0.1 mg/ml.

Description

Prazobind is an α_1 -adrenergic receptor (AR) antagonist.¹ It inhibits binding of the α_1 -AR ligand, BE-2254, in rat hippocampus and liver, which highly express α_{1A} - and α_{1B} -ARs, respectively (IC₅₀ = 1 nM for both).² It decreases norepinephrine-induced contraction of isolated guinea pig aorta when used at a concentration of 1 nM.³

References

1. Helman, J., Kusiak, J.W., Pitha, J., *et al.* Inhibition of α_1 -adrenergic responsiveness in intact cells by a new, irreversible receptor antagonist. *Biochem. Biophys. Res. Commun.* **142**(2), 403-409 (1987).
2. Mante, S. and Minneman, K.P. The alkylating prazosin analog SZL 49 inactivates both α_{1A} - and α_{1B} -adrenoceptors. *Eur. J. Pharmacol.* **208**(2), 113-117 (1991).
3. Oriowo, M. α_1 -adrenoceptor subtype(s) mediating noradrenaline-induced contractions of the guinea-pig aorta. *Fundam. Clin. Pharmacol.* **8**(3), 214-219 (1994).

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, 10/17/2022

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