

PRODUCT INFORMATION



CGS 9343B

Item No. 32844

CAS Registry No.: 109826-27-9
Formal Name: 1,3-dihydro-1-[1-[(4-methyl-4H,6H-pyrrolo[1,2-a][4,1]benzoxazepin-4-yl)methyl]-4-piperidinyl]-2H-benzimidazol-2-one, 2Z-butenedioate

Synonyms: KW-5617, Zaldaride Maleate

MF: C₂₆H₂₈N₄O₂ • C₄H₄O₄

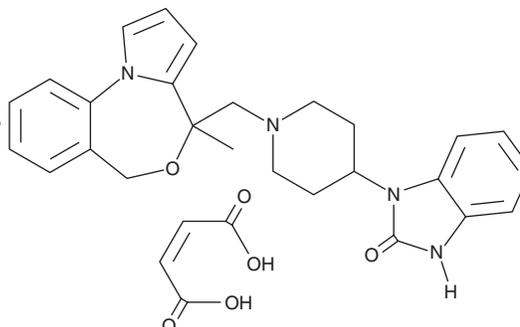
FW: 544.6

Purity: ≥98%

Supplied as: A 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

CGS 9343B is supplied as a solid. A stock solution may be made by dissolving the CGS 9343B in the solvent of choice, which should be purged with an inert gas. CGS 9343B is soluble in DMSO and slightly soluble in ethanol (warmed).

Description

CGS 9343B is a calmodulin inhibitor.¹⁻³ It inhibits calmodulin-induced phosphodiesterase (PDE) activity (IC₅₀ = 3.3 μM) but does not inhibit basal PKC activity at 100 μM.² CGS 9343B (120 μM) inhibits calcium release induced by inositol 1,4,5-trisphosphate (IP₃) in 261B rat liver epithelial cells.² It reduces the incidence of diarrhea without affecting gastrointestinal transit time in a mouse model of secretory diarrhea induced by castor oil when administered at a dose of 3 mg/kg.³

References

1. Norman, J.A., Ansell, J., Stone, G.A., *et al.* CGS 9343B, a novel, potent, and selective inhibitor of calmodulin activity. *Mol. Pharmacol.* **31(5)**, 535-540 (1987).
2. Hill, T.D., Campos-Gonzalez, R., Kindmark, H., *et al.* Inhibition of inositol trisphosphate-stimulated calcium mobilization by calmodulin antagonists in rat liver epithelial cells. *J. Biol. Chem.* **263(31)**, 16479-16484 (1988).
3. Shook, J.E., Burks, T.F., Wasley, J.W., *et al.* Novel calmodulin antagonist CGS 9343B inhibits secretory diarrhea. *J. Pharmacol. Exp. Ther.* **251(1)**, 247-252 (1989).

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, 11/08/2022

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