

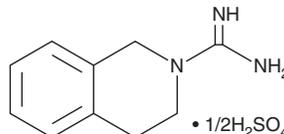
PRODUCT INFORMATION



Debrisoquin (hemisulfate)

Item No. 29568

CAS Registry No.: 581-88-4
Formal Name: 3,4-dihydro-2(1H)-isoquinolinecarboximidamide, hemisulfate
Synonyms: Isocaramidine sulfate, Ro 5-3307/1
MF: C₁₀H₁₃N₃ • 1/2H₂SO₄
FW: 224.3
Purity: ≥98%
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

Debrisoquin (hemisulfate) is supplied as a crystalline solid. Aqueous solutions of debrisoquin (hemisulfate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of debrisoquin (hemisulfate) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Debrisoquin is a post-ganglionic sympathetic blocker.¹ It increases accumulation of the monoamine oxidase (MAO) substrate *l*-*m*-octopamine in isolated rabbit heart when used at a concentration of 1 mM.² Debrisoquin reduces norepinephrine levels in the left and right atrium and ventricles, as well as the mesenteric and femoral arteries in anesthetized dogs when administered at a dose of 5 mg/kg per day for seven days.³ Debrisoquin decreases mean aortic pressure, heart rate, and total peripheral vascular resistance in the same model.

References

1. Heffernan, A.G.A. and Carty, A.T. Clinical observations on the use of debrisoquine sulphate (declinax) in the treatment of hypertension. *I. J. Med. Sc.* **3(1)**, 37-43 (1970).
2. Giachetti, A. and Shore, P.A. Monoamine oxidase inhibition in the adrenergic neuron by bretylium, debrisoquin, and other adrenergic neuronal blocking agents. *Biochem. Pharmacol.* **16(1)**, 237-238 (1967).
3. Cavero, I., Gerold, M., Saner, A., et al. The cardiovascular effects of the antihypertensive drug debrisoquin: A contribution to the pharmacology of chronic treatment. I. One-week administration to dogs. *J. Pharmacol. Exp. Ther.* **206(1)**, 123-131 (1978).

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, 09/30/2022

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