

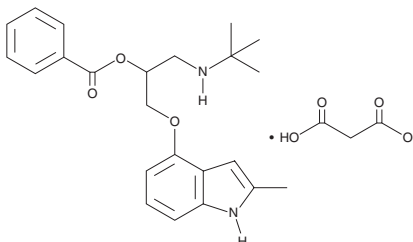
PRODUCT INFORMATION



Bopindolol (malonate)

Item No. 23365

CAS Registry No.: 82857-38-3
Formal Name: propanedioic acid compound with 2-[(1,1-dimethylethyl)amino]-1-[[[(2-methyl-1H-indol-4-yl)oxy)methyl]ethyl]benzoate
MF: C₂₃H₂₈N₂O₃ • C₃H₄O₄
FW: 484.5
Purity: ≥98%
Supplied as: A solid
Storage: Room temperature
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Bopindolol (malonate) is supplied as a solid. A stock solution may be made by dissolving the bopindolol (malonate) in water. The solubility of bopindolol (malonate) in water is approximately 5 mM. We do not recommend storing the aqueous solution for more than one day.

Description

Bopindolol is a non-selective β -adrenergic receptor antagonist.¹ *In vivo*, bopindolol decreases heart rate and blood pressure and attenuates ischemia-induced myocardial acidosis in dogs.² Formulations containing bopindolol have been used to treat essential and renovascular hypertension.¹

References

1. van Brummelen, P., Bühler, F.R., Amann, F.W., *et al.* Effects of a new long-acting beta-blocker bopindolol (LT 31-200) on blood pressure, plasma catecholamines, renin and cholesterol in patients with arterial hypertension. *Eur. J. Pharmacol.* **22(6)**, 491-493 (1982).
2. Abiko, Y., Gotoh, H., Yokoyama, T., *et al.* Bopindolol and its metabolite 18-502 attenuate regional myocardial acidosis during partial occlusion of the coronary artery in dogs. *Arch. Int. Pharmacodyn. Ther.* **327(1)**, 40-45 (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