

PRODUCT INFORMATION

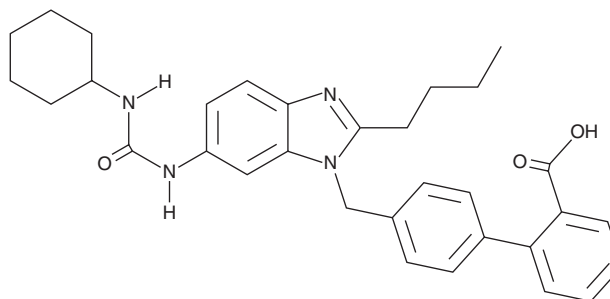


BIBS39

Item No. 22932

CAS Registry No.: 133085-33-3
Formal Name: 4'-[[2-butyl-6-[[cyclohexylamino]carbonyl]amino]-1H-benzimidazol-1-yl]methyl]-[1,1'-biphenyl]-2-carboxylic acid

MF: C₃₂H₃₆N₄O₃
FW: 524.7
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 295 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

BIBS39 is supplied as a crystalline solid. A stock solution may be made by dissolving the BIBS39 in the solvent of choice, which should be purged with an inert gas. BIBS39 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BIBS39 in these solvents is approximately 30 mg/ml.

BIBS39 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BIBS39 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BIBS39 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BIBS39 is a nonpeptide antagonist of the angiotensin II type 1 (AT₁) receptor (K_i = 29 nM in rat lung membranes).¹ It is selective for AT₁ over AT₂ receptors (K_i = 480 nM in isolated rat adrenal medulla), as well as α₁-, α₂-, β₁-, and β₂-adrenergic, adenosine A₁ and A₂, muscarinic M₁, M₂, and M₃, and atrial natriuretic peptide receptors (K_is = >100 μM for all). BIBS39 (0.3-10 mg/kg) decreases diastolic blood pressure in pithed normotensive rats without affecting heart rate and in renal hypertensive rats with an increase in heart rate when administered at doses ranging from 1 to 30 mg/kg.^{1,2}

References

1. Zhang, J., Entzeroth, M., Wiene, W., *et al.* Characterization of BIBS 39 and BIBS 222: Two new nonpeptide angiotensin II receptor antagonists. *Eur. J. Pharmacol.* **218(1)**, 35-41 (1992).
2. Zhang, J., Van Meel, J.C., and Van Zwieten, P.A. Antihypertensive activity of the novel, nonpeptide angiotensin II receptor antagonists BIBS 39 and BIBS 222 in conscious renal hypertensive rats. *Pharmacology* **46(5)**, 241-247 (1993).

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, 11/14/2022

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