

PRODUCT INFORMATION



BIBO-3304 (trifluoroacetate salt)

Item No. 35205

CAS Registry No.: 191868-14-1
Formal Name: N-[(1R)-1-[[[4-[[[aminocarbonyl]amino]methyl]phenyl]methyl]amino]carbonyl]-4-[(aminoiminomethyl)amino]butyl]- α -phenyl-benzeneacetamide, trifluoroacetate salt

MF: C₂₉H₃₅N₇O₃ • CF₃COOH

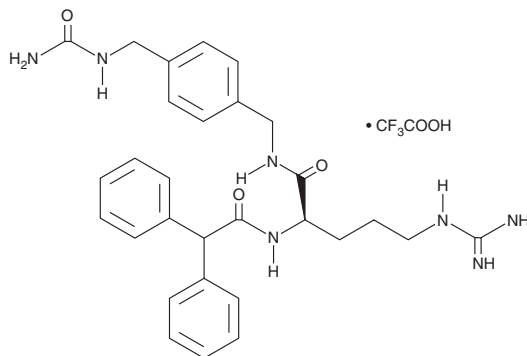
FW: 643.7

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

BIBO-3304 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the BIBO-3304 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. BIBO-3304 (trifluoroacetate salt) is soluble in the organic solvent DMSO.

Description

BIBO-3304 is an antagonist of the neuropeptide Y (NPY) receptor Y₁ (K_is = 0.38 and 0.72 nM, respectively).¹ It is selective for Y₁ over Y₂, Y₄, and Y₅ receptors (K_is = >1,000 nM for all). BIBO-3304 (30 µg/animal) inhibits NPY-induced increases in food intake in fasted rats when injected into the paraventricular nucleus (PVN) of the hypothalamus. It also inhibits the pressor response induced by NPY or bilateral common carotid artery occlusion in cats when administered at a dose of 100 µg/kg.²

References

1. Wieland, H.A., Engel, W., Eberlein, W., *et al.* Subtype selectivity of the novel nonpeptide neuropeptide Y Y₁ receptor antagonist BIBO 3304 and its effect on feeding in rodents. *Br. J. Pharmacol.* **125**(3), 549-555 (1998).
2. Capurro, D. and Huidobro-Toro, J.P. The involvement of neuropeptide Y Y₁ receptors in the blood pressure baroreflex: Studies with BIBP 3226 and BIBO 3304. *Eur. J. Pharmacol.* **376**(3), 251-255 (1999).

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, 10/11/2022

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