

PRODUCT INFORMATION

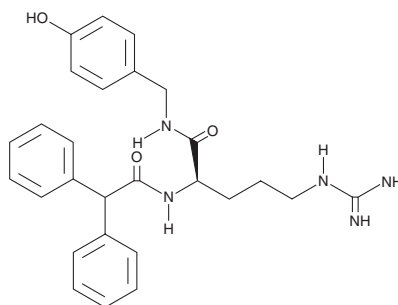


BIBP3226

Item No. 33210

CAS Registry No.: 159013-54-4
Formal Name: N-[(1R)-4-[(aminoiminomethyl)amino]-1-[[[4-hydroxyphenyl)methyl]amino]carbonyl]butyl]- α -phenylbenzeneacetamide

MF: C₂₇H₃₁N₅O₃
FW: 473.6
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

BIBP3226 is supplied as a solid. A stock solution may be made by dissolving the BIBP3226 in the solvent of choice, which should be purged with an inert gas. BIBP3226 is soluble in ethanol and DMSO.

Description

BIBP3226 is a nonpeptide antagonist of neuropeptide Y (NPY) receptor Y₁ (K_i = 1.1 nM).¹ It is selective for Y₁ over Y₂, Y₄, and Y₅ receptors (K_is = >1,000 nM for all). It also binds to neuropeptide FF receptor 1 (NPFF1) and NPFF2 (K_is = 108 and 79 nM, respectively) and reverses NPFF-induced inhibition of forskolin-induced cAMP accumulation in CHO cells expressing human NPFF2 in a concentration-dependent manner.¹ BIBP3226 inhibits NPY-induced increases in perfusion pressure in isolated rat kidney but not the NPY-induced twitch response in isolated rat vas deferens (IC₅₀s = 26 and >10,000 nM, respectively).² BIBP3226 inhibits NPY-induced increases in blood pressure in pithed rats (ED₅₀ = 0.11 mg/kg).² It also inhibits NPFF-induced hypothermia in mice when administered intracerebroventricularly (i.c.v.) at a dose of 5 nmol.³

References

1. Mollereau, C., Gouardères, C., Dumont, Y., *et al.* Agonist and antagonist activities on human NPFF₂ receptors of the NPY ligands GR231118 and BIBP3226. *Br. J. Pharmacol.* **133**(1), 1-4 (2001).
2. Rudolf, K., Eberlein, W., Engel, W., *et al.* The first highly potent and selective non-peptide neuropeptide Y Y₁ receptor antagonist: BIBP3226. *Eur. J. Pharmacol.* **271**(2-3), R11-R13 (1994).
3. Fang, Q., Guo, J., He, F., *et al.* In vivo inhibition of neuropeptide FF agonism by BIBP3226, an NPY Y1 receptor antagonist. *Peptides* **27**(9), 2207-2213 (2006).

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, 12/05/2022

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