

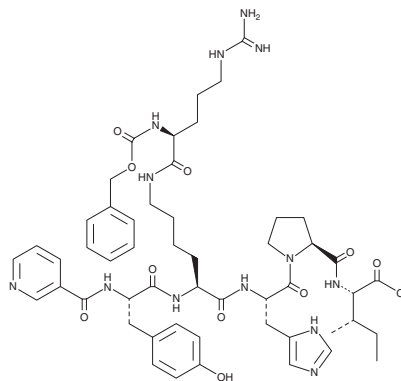
PRODUCT INFORMATION



CGP 42112

Item No. 32874

CAS Registry No.: 127060-75-7
Formal Name: N-(3-pyridinylcarbonyl)-L-tyrosyl-N⁶-[N²-[(phenylmethoxy)carbonyl]-L-arginyl]-L-lysyl-L-histidyl-L-prolyl-L-isoleucine
MF: C₅₂H₆₉N₁₃O₁₁
FW: 1,052.2
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

CGP 42112 is supplied as a crystalline solid. A stock solution may be made by dissolving the CGP 42112 in the solvent of choice, which should be purged with an inert gas. CGP 42112 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CGP 42112 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CGP 42112 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of CGP 42112 in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

CGP 42112 is an angiotensin II type 2 (AT₂) receptor agonist.^{1,2} It binds to AT₂ receptors (K_i = 0.24 nM) and inhibits cGMP production and tyrosine hydroxylase activity in primary porcine adrenal medullary cells when used at concentrations ranging from 0.001 to 1 μM, effects that can be blocked by the AT₂ receptor antagonist PD 123319 (Item No. 16099).

References

1. Criscione, L., Thomann, H., Whitebread, S., *et al.* Binding characteristics and vascular effects of various angiotensin II antagonists. *J. Cardiovasc. Pharmacol.* **16 (Suppl 4)**, S56-S59 (1990).
2. Takekoshi, K., Isobe, K.I., Nanmoku, T., *et al.* Angiotensin-II subtype 2 receptor agonist (CGP-42112) inhibits catecholamine biosynthesis in cultured porcine adrenal medullary chromaffin cells. *Biochem. Biophys. Res. Commun.* **272(2)**, 544-550 (2000).

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, 11/11/2022

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