

PRODUCT INFORMATION



CGP 20712A

Item No. 40765

CAS Registry No.: 105737-62-0

Formal Name: 2-hydroxy-5-[2-[[2-hydroxy-3-[4-[1-methyl-4-(trifluoromethyl)-1H-imidazol-2-yl]phenoxy]propyl]amino]ethoxy]-benzamide, monomethanesulfonate

MF: $C_{23}H_{25}F_3N_4O_5 \cdot CH_3SO_3H$

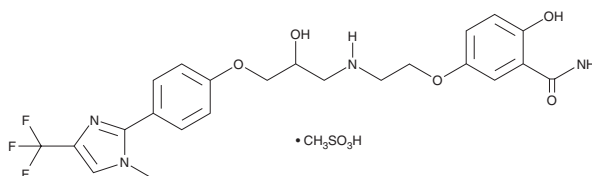
FW: 590.6

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

CGP 20712A is supplied as a solid. A stock solution may be made by dissolving the CGP 20712A in the solvent of choice, which should be purged with an inert gas. CGP 20712A is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in methanol and acetonitrile.

Description

CGP 20712A is an antagonist of the β_1 -adrenergic receptor (β_1 -AR; $IC_{50} = 0.7$ nM).¹ It is selective for β_1 -ARs over β_2 -ARs ($IC_{50} = 6,700$ nM). CGP 20712 A (0.1-1,000 μM) inhibits adenylyl cyclase activity induced by the β_1 - and β_2 -AR agonist isoproterenol (Item No. 15592) in brown adipose tissue plasma membranes isolated from both lean and pre-obese rat pups.²

References

1. Dooley, D.J., Bittiger, H., and Reymann, N.C. CGP 20712 A: A useful tool for quantitating β_1 - and β_2 -adrenoceptors. *Eur. J. Pharmacol.* **130(1-2)**, 137-139 (1986).
2. Charon, C., Krief, S., Diot-Dupuy, F., et al. Early alterations in the brown adipose tissue adenylyl cyclase system of pre-obese Zucker rat *fa/fa* pups: Decreased G-proteins and β_3 -adrenoceptor activities. *Biochem J.* **312(Pt 3)**, 781-788 (1995)

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, 04/01/2024

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