

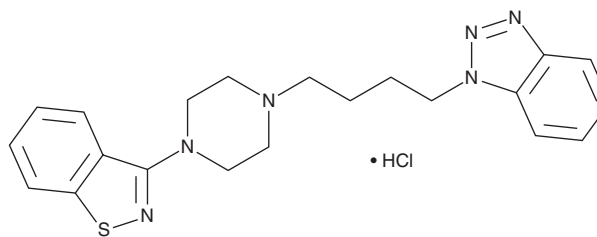
PRODUCT INFORMATION



MT-1207 (hydrochloride)

Item No. 41720

CAS Registry No.: 1610794-70-1
Formal Name: 1-[4-[4-(1,2-benzisothiazol-3-yl)-1-piperazinyl]butyl]-1H-benzotriazole, monohydrochloride
MF: $C_{21}H_{24}N_6S \cdot HCl$
FW: 429.0
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

MT-1207 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the MT-1207 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. MT-1207 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

MT-1207 is an antagonist of α_{1A} -, α_{1B} -, and α_{1D} -adrenergic receptors and the serotonin (5-HT) receptor subtype 5-HT_{2A} (IC_{50} s = <0.1, 0.15, 1.4, and 0.27 nM, respectively).¹ It is selective for these receptors over a panel of 87 receptors, ion channels, and enzymes at 1 μ M but does inhibit α_{2A} - and α_{2B} -adrenergic and 5-HT_{1A}, 5-HT_{1B}, 5-HT_{2B}, and 5-HT_{2C} receptors (IC_{50} s = 18.1, 4.89, 4.68, 41, 11.5, and 26 nM, respectively), among others. MT-1207 induces relaxation of isolated rat aortic rings precontracted with epinephrine, potassium chloride, norepinephrine, or 5-HT in a concentration-dependent manner. It decreases systolic and diastolic blood pressure and heart rate in spontaneously hypertensive rats and two kidney-one clip (2K-1C) renal hypertensive dogs when administered at doses of 10 and 4 mg/kg, respectively.

Reference

1. Xu, T.Y., Wang, P., Tian, J.S., *et al.* Pharmacological characterization of MT-1207, a novel multitarget antihypertensive agent. *Acta. Pharmacol. Sin.* **42(6)**, 885-897 (2021).

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, 10/14/2024

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