

PRODUCT INFORMATION



Atrasentan (hydrochloride)

Item No. 31238

CAS Registry No.: 195733-43-8

Formal Name: 4-(1,3-benzodioxol-5-yl)-1-[2R-(dibutylamino)-2-oxoethyl]-2-(4S-methoxyphenyl)-3R-pyrrolidinecarboxylic acid, monohydrochloride
A-127722, A-147627, ABT-627

Synonyms:

MF: $C_{29}H_{38}N_2O_6 \cdot HCl$

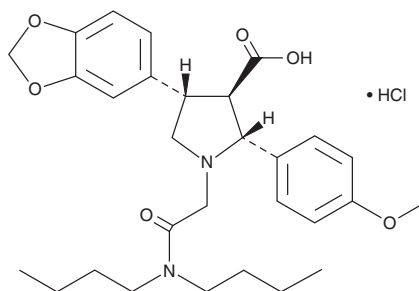
FW: 547.1

Purity: $\geq 98\%$

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

Atrasentan (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the atrasentan (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Atrasentan (hydrochloride) is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Atrasentan is an antagonist of endothelin receptor type A (ET_A ; $K_i = 69$ pM for the human receptor).¹ It is selective for ET_A over ET_B ($K_i = 139$ nM for the human receptor) and a panel of 29 other receptors at 10 μM . Atrasentan inhibits endothelin-1-induced secretion of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) from human pericardial smooth muscle cells (HPSMCs; $pA_2 = 10.5$), as well as endothelin-1-induced vasoconstriction in isolated rat aortic rings ($pA_2 = 9.2$). It inhibits endothelin-1-induced increases in the pressor response in normotensive rats when administered at a dose of 10 mg/kg. Atrasentan (5 mg/kg per day) reduces urinary levels of albumin, TGF- β , and a prostaglandin E_2 metabolite, as well as decreases the number of macrophages in the renal cortex, in a rat model of streptozotocin-induced diabetic nephropathy.²

References

1. Opgenorth, T.J., Adler, A.L., Calzadilla, S.V., *et al.* Pharmacological characterization of A-127722: An orally active and highly potent ET_A -selective receptor antagonist. *J. Pharmacol. Exp. Ther.* **276**(2), 473-481 (1996).
2. Sasser, J.M., Sullivan, J.C., Hobbs, J.L., *et al.* Endothelin A receptor blockade reduces diabetic renal injury via an anti-inflammatory mechanism. *J. Am. Soc. Nephrol.* **18**(1), 143-154 (2007).

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, 08/07/2024

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