

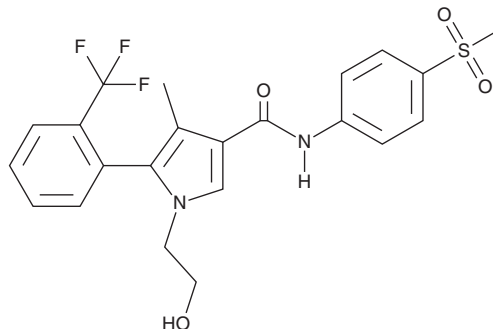
PRODUCT INFORMATION



Esaxerenone

Item No. 41082

CAS Registry No.: 1632006-28-0
Formal Name: 1-(2-hydroxyethyl)-4-methyl-N-[4-(methylsulfonyl)phenyl]-5S-[2-(trifluoromethyl)phenyl]-1H-pyrrole-3-carboxamide
Synonym: CS-3150
MF: C₂₂H₂₁F₃N₂O₄S
FW: 466.5
Purity: ≥98%
Supplied as: A 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

Esaxerenone is supplied as a solid. A stock solution may be made by dissolving the esaxerenone in the solvent of choice, which should be purged with an inert gas. Esaxerenone is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

Esaxerenone is a non-steroidal mineralocorticoid receptor antagonist (IC₅₀ = 9.4 nM).¹ It is selective for mineralocorticoid receptors over glucocorticoid, androgen, and progesterone receptors (IC₅₀s = >10,000 nM for all). Esaxerenone (0.3, 1, or 3 mg/kg) increases the urinary ratio of sodium to potassium in bilateral adrenalectomized rats. It decreases systolic blood pressure in DOCA-salt hypertensive rats when administered at a dose of 3 mg/kg to a greater extent than the mineralocorticoid receptor antagonists spironolactone or eplerenone (Item No. 15616). Esaxerenone (2 mg/kg per day) reduces systolic blood pressure, urinary protein levels, blood levels of brain natriuretic peptide (BNP), kidney and left ventricle weights, and glomerulosclerosis, as well as increases blood potassium levels, in Dahl salt-sensitive rats fed a high-salt diet.² Formulations containing esaxerenone have been used in the treatment of hypertension.

References

1. Arai, K., Homma, T., Morikawa, Y., *et al.* Pharmacological profile of CS-3150, a novel, highly potent and selective non-steroidal mineralocorticoid receptor antagonist. *Eur. J. Pharmacol.* **761**, 226-234 (2015).
2. Arai, K., Tsuruoka, H., and Homma, T. CS-3150, a novel non-steroidal mineralocorticoid receptor antagonist, prevents hypertension and cardiorenal injury in Dahl salt-sensitive hypertensive rats. *Eur. J. Pharmacol.* **769**, 266-273 (2015).

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, 07/19/2024

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