

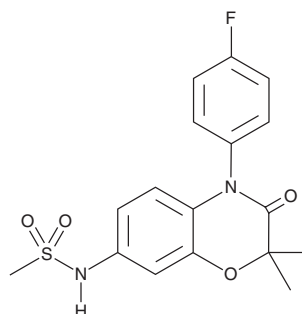
PRODUCT INFORMATION



Apararenone

Item No. 30671

CAS Registry No.: 945966-46-1
Formal Name: N-[4-(4-fluorophenyl)-3,4-dihydro-2,2-dimethyl-3-oxo-2H-1,4-benzoxazin-7-yl]-methanesulfonamide
Synonym: MT-3995
MF: C₁₇H₁₇FN₂O₄S
FW: 364.4
Purity: ≥95%
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

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

Description

Apararenone is a non-steroidal mineralocorticoid receptor antagonist ($K_i = 0.104 \mu\text{M}$ for the human receptor).¹ It is selective for mineralocorticoid receptors over estrogen receptor α (ER α) and ER β and glucocorticoid, progesterone, and androgen receptors ($\text{IC}_{50}\text{s} = >100 \mu\text{M}$ for all). Apararenone increases the urinary ratio of sodium to potassium in adrenalectomized rats in a dose-dependent manner. It inhibits aldosterone- and nephrectomy-induced increases in systolic blood pressure and urinary albumin levels, as well as myocardial necrosis and fibrosis, cardiac arterial degeneration, and renal tubule damage, in a rat model of renal hypertension when administered at doses of 3 or 10 mg/kg per day.

Reference

1. Iijima, T., Katoh, M., Takedomi, K., et al. Discovery of apararenone (MT-3995) as a highly selective, potent, and novel nonsteroidal mineralocorticoid receptor antagonist. *J. Med. Chem.* **65**(12), 8127-8143 (2022).

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, 10/23/2024

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