

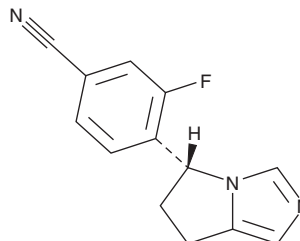
PRODUCT INFORMATION



Osilodrostat

Item No. 33928

CAS Registry No.: 928134-65-0
Formal Name: 4-[(5R)-6,7-dihydro-5H-pyrrolo[1,2-c]imidazol-5-yl]-3-fluoro-benzonitrile
Synonym: LCI699
MF: $C_{13}H_{10}FN_3$
FW: 227.2
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 227 nm
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

Osilodrostat is supplied as a solid. A stock solution may be made by dissolving the osilodrostat in the solvent of choice, which should be purged with an inert gas. Osilodrostat is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of osilodrostat in these solvents is approximately 30 mg/ml.

Description

Osilodrostat is an inhibitor of aldosterone synthase ($IC_{50} = 0.7$ nM for the recombinant human enzyme).¹ It is selective for human aldosterone synthase over rat aldosterone synthase ($IC_{50} = 160$ nM). Osilodrostat inhibits increases in plasma aldosterone levels induced by angiotensin II or adrenocorticotrophic hormone (ACTH) in rats (ED_{50} s = 0.6 and 1.1 mg/kg, respectively). It reduces mean arterial pressure and prolongs survival in rats expressing human renin and angiotensinogen, a double-transgenic model of angiotensin II-dependent hypertension, in a dose-dependent manner. Formulations containing osilodrostat have been used in the treatment of Cushing's syndrome.

Reference

1. Ménard, J., Rigel, D.F., Watson, C., *et al.* Aldosterone synthase inhibition: Cardiorenal protection in animal disease models and translation of hormonal effects to human subjects. *J. Transl. Med.* **12**, 340 (2014).

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, 12/01/2022

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