

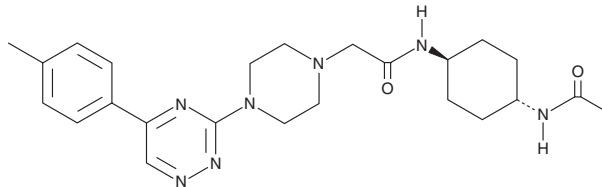
PRODUCT INFORMATION



Lorundrostat

Item No. 43548

CAS Registry No.: 1820940-17-7
Formal Name: N-[*trans*-4-(acetylamino)cyclohexyl]-4-[5-(4-methylphenyl)-1,2,4-triazin-3-yl]-1-piperazineacetamide
Synonyms: MLS-101, MT-4129
MF: C₂₄H₃₃N₇O₂
FW: 451.6
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

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

Description

Lorundrostat is an inhibitor of aldosterone synthase, also known as cytochrome P450 (CYP) 11B2 (CYP11B2; IC₅₀ = 1.27 nM).¹ It is selective for aldosterone synthase over CYP11B1 (IC₅₀ = 475 nM).

Reference

1. Shimizu, H., Tortorici, M.A., Ohta, Y., *et al.* First-in-human study evaluating safety, pharmacokinetics, and pharmacodynamics of lorundrostat, a novel and highly selective aldosterone synthase inhibitor. *Clin Transl Sci.* **17**(8), e70000 (2024).

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, 07/22/2025

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