

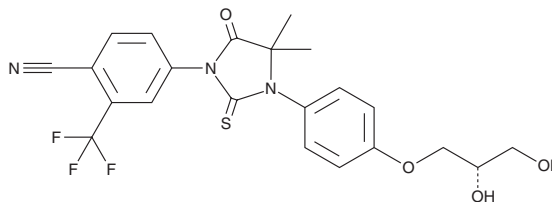
PRODUCT INFORMATION



Rezvilutamide

Item No. 45332

CAS Registry No.: 1572045-62-5
Formal Name: 4-[3-[4-[(2S)-2,3-dihydroxypropoxy]phenyl]-4,4-dimethyl-5-oxo-2-thioxo-1-imidazolidinyl]-2-(trifluoromethyl)-benzonitrile
Synonym: SHR-3680
MF: C₂₂H₂₀F₃N₃O₄S
FW: 479.5
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

Rezvilutamide is supplied as a solid. A stock solution may be made by dissolving the rezvilutamide in the solvent of choice, which should be purged with an inert gas. Rezvilutamide is soluble (≥10 mg/ml) in DMSO. Rezvilutamide is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Rezvilutamide is a non-steroidal androgen receptor antagonist.¹ It reduces secretion of prostate-specific androgen (PSA) from LNCaP prostate cancer cells (IC₅₀ = 22 nM) and reduces proliferation of the same cells (IC₅₀ = 33 nM). Formulations containing rezvilutamide have been used in the treatment of prostate cancer.

Reference

1. Lu, H., Sun, P., Fei, H., *et al.* Imidazoline derivatives, preparation methods thereof, and their applications in medicine. *Shanghai Hengrui Pharmaceutical Co., LTD., WO2014036897* (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, 04/17/2026

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