

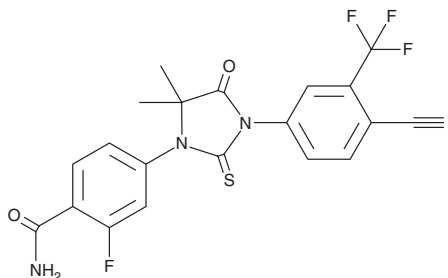
PRODUCT INFORMATION



N-desmethyl Enzalutamide

Item No. 41025

CAS Registry No.: 1242137-16-1
Formal Name: 4-[3-[4-cyano-3-(trifluoromethyl)phenyl]-5,5-dimethyl-4-oxo-2-thioxo-1-imidazolidinyl]-2-fluoro-benzamide
Synonym: N-desmethyl MDV 3100
MF: C₂₀H₁₄F₄N₄O₂S
FW: 450.4
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

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

Description

N-desmethyl Enzalutamide is an active metabolite of the androgen receptor antagonist enzalutamide (Item No. 11596).¹ It is formed from enzalutamide by the cytochrome P450 (CYP) isoforms CYP3A4 and CYP2C8.²

References

1. Gibbons, J.A., Ouatas, T., Krauwinkel, W., *et al.* Clinical pharmacokinetic studies of enzalutamide. *Clin. Pharmacokinet.* **54**(10), 1043-1055 (2015).
2. Gibbons, J.A., de Vries, M., Krauwinkel, W., *et al.* Pharmacokinetic drug interaction studies with enzalutamide. *Clin. Pharmacokinet.* **54**(10), 1057-1069 (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, 07/10/2024

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