

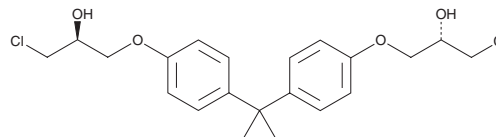
PRODUCT INFORMATION



Ralaniten

Item No. 42391

CAS Registry No.: 1203490-23-6
Formal Name: (2R)-3-[4-[1-[4-[(2S)-3-chloro-2-hydroxypropoxy]phenyl]-1-methylethyl]phenoxy]-1,2-propanediol
Synonym: EPI-002
MF: C₂₁H₂₇ClO₅
FW: 394.9
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

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

Description

Ralaniten is an antagonist of the androgen receptor N-terminal domain (NTD).¹ It inhibits androgen receptor NTD transactivation in a reporter assay (IC₅₀ = 7.4 μM). Ralaniten (10-30 μM) reduces androgen-dependent growth of LNCaP prostate cancer cells. *In vivo*, ralaniten (50 mg/kg) decreases tumor volume in an LNCaP castration-resistant prostate cancer (CRPC) mouse xenograft model.

Reference

1. Myung, J.K., Banuelos, C.A., Fernandez, J.G., *et al.* An androgen receptor N-terminal domain antagonist for treating prostate cancer. *J. Clin. Invest.* **123**(7), 2948-2960 (2013).

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, 12/17/2024

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