

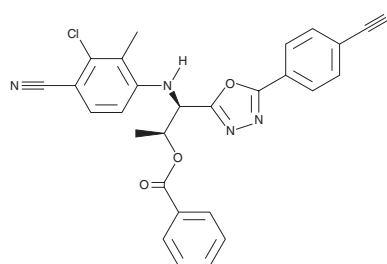
PRODUCT INFORMATION



RAD150

Item No. 39173

CAS Registry No.: 1208070-53-4
Formal Name: 4-[[[(1R,2S)-2-(benzoyloxy)-1-[5-(4-cyanophenyl)-1,3,4-oxadiazol-2-yl]propyl]amino]-2-chloro-3-methyl-benzonitrile
MF: C₂₇H₂₀ClN₅O₃
FW: 497.9
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

RAD150 is supplied as a solid. A stock solution may be made by dissolving the RAD150 in the solvent of choice, which should be purged with an inert gas. RAD150 is soluble in methanol.

Description

RAD150 is a nonsteroidal androgen receptor modulator (IC₅₀ = 130 nM) and benzoate ester derivative of the selective androgen receptor modulator (SARM) RAD140 (Item No. 18773).¹

Reference

1. Miller, C.P. Selective androgen receptor modulators. *Radius Health, Inc. US2010/0041721A1* (2010).

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, 01/02/2024

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