

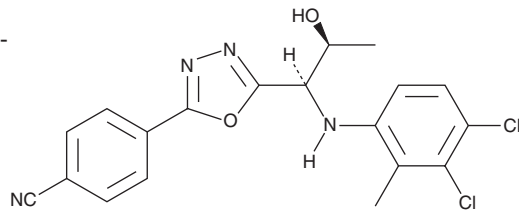
PRODUCT INFORMATION



RAD140

Item No. 18773

CAS Registry No.: 1182367-47-0
Formal Name: 2-chloro-4-[[[(1R,2S)-1-[5-(4-cyanophenyl)-1,3,4-oxadiazol-2-yl]-2-hydroxypropyl]amino]-3-methyl-benzonitrile
Synonym: Testolone
MF: C₂₀H₁₆ClN₅O₂
FW: 393.8
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 277 nm
Supplied as: A crystalline 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

RAD140 is supplied as a crystalline solid. A stock solution may be made by dissolving the RAD140 in the solvent of choice, which should be purged with an inert gas. RAD140 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of RAD140 in ethanol is approximately 2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

RAD140 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RAD140 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. RAD140 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

RAD140 is an orally bioavailable, nonsteroidal selective androgen receptor modulator (SARM) that binds to the androgen receptor with a K_i value of 7 nM.¹ At 0.03-0.3 mg/kg, it has been shown to stimulate muscle weight increase without affecting prostate or seminal vesicle in castrated male rats.¹ At 0.3-1 mg/kg, RAD140 demonstrates partial antagonist activity in seminal vesicle and prostate.¹

Reference

1. Basaria, S., Collins, L., Dillon, E.L., *et al.* The safety, pharmacokinetics, and effects of LGD-4033, a novel nonsteroidal oral, selective androgen receptor modulator, in healthy young men. *J. Gerontol. A Biol. Sci. Med. Sci.* **68(1)**, 87-95 (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, 11/07/2022

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