

PRODUCT INFORMATION

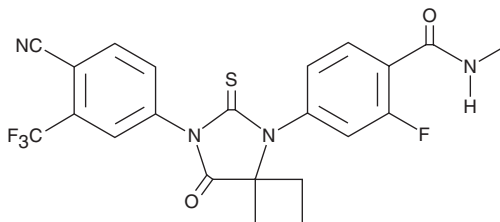


RD162

Item No. 13039

CAS Registry No.: 915087-27-3
Formal Name: 4-[7-[4-cyano-3-(trifluoromethyl)phenyl]-8-oxo-6-thioxo-5,7-diazaspiro[3.4]oct-5-yl]-2-fluoro-N-methyl-benzamide

MF: C₂₂H₁₆F₄N₄O₂S
FW: 476.4
Purity: ≥98%
UV/Vis.: λ_{max}: 239 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

RD162 is supplied as a crystalline solid. A stock solution may be made by dissolving the RD162 in the solvent of choice, which should be purged with an inert gas. RD162 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of RD162 in these solvents is approximately 30 mg/ml. RD162 is slightly soluble in ethanol.

RD162 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RD162 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. RD162 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

RD162 is a second-generation androgen receptor (AR) antagonist. It inhibits binding of ¹⁸F-FDHT to AR in human prostate cancer cells with an IC₅₀ value of 30.9 nM.¹ RD162 completely suppresses VCaP cell proliferation at 1 μM and induces apoptosis of VCaP cells at 10 μM. RD162 is bioavailable and reduces LNCaP/AR tumor growth in castrated male mice at a daily oral dose of 10 mg/kg.¹

Reference

1. Tran, C., Ouk, S., Clegg, N.J., *et al.* Development of a second-generation antiandrogen for treatment of advanced prostate cancer. *Science* **324**(5928), 787-790 (2009).

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, 11/02/2022

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