

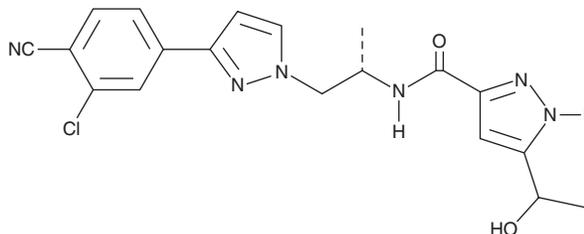
PRODUCT INFORMATION



ODM-201

Item No. 25643

CAS Registry No.: 1297538-32-9
Formal Name: N-[(1S)-2-[3-(3-chloro-4-cyanophenyl)-1H-pyrazol-1-yl]-1-methylethyl]-5-(1-hydroxyethyl)-1H-pyrazole-3-carboxamide
Synonym: BAY 1841788
MF: C₁₉H₁₉ClN₆O₂
FW: 398.9
Purity: ≥98%
UV/Vis.: λ_{max}: 288 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

ODM-201 is supplied as a crystalline solid. A stock solution may be made by dissolving the ODM-201 in the solvent of choice, which should be purged with an inert gas. ODM-201 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ODM-201 in ethanol is approximately 10 mg/ml and approximately 25 mg/ml in DMSO and DMF.

ODM-201 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ODM-201 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ODM-201 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ODM-201 is an androgen receptor (AR) antagonist ($K_i = 11$ nM).¹ It inhibits AR transactivation in U2OS osteosarcoma cells expressing human wild-type and mutant ARs (IC_{50} s = 65, 66, 1,500, and 1,782 nM for wild-type, AR^{F876L}, AR^{W741L}, and AR^{T877A}, respectively). ODM-201 prevents androgen-induced AR nuclear translocation in AR-overexpressing HS-HEK293 and LNCaP cells and suppresses androgen-induced proliferation of VCaP cells ($IC_{50} = 230$ nM). *In vivo*, ODM-201 (50 mg/kg per day) reduces tumor growth in a VCaP castrated mouse xenograft model. It also inhibits tumor growth without increasing serum testosterone levels in a VCaP intact mouse xenograft model.

Reference

1. Moilanen, A.M., Riikonen, R., Oksala, R., *et al.* Discovery of ODM-201, a new-generation androgen receptor inhibitor targeting resistance mechanisms to androgen signaling-directed prostate cancer therapies. *Sci. Rep.* 5:12007 (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, 12/07/2022

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