

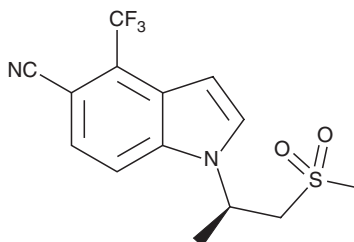
PRODUCT INFORMATION



GSK2881078

Item No. 20313

CAS Registry No.: 1539314-06-1
Formal Name: 1-[(1R)-1-methyl-2-(methylsulfonyl)ethyl]-4-(trifluoromethyl)-1H-indole-5-carbonitrile
MF: $C_{14}H_{13}F_3N_2O_2S$
FW: 330.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 243, 303 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GSK2881078 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2881078 in the solvent of choice, which should be purged with an inert gas. GSK2881078 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GSK2881078 in ethanol is approximately 20 mg/ml and approximately 25 mg/ml in DMSO and DMF.

GSK2881078 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK2881078 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK2881078 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK2881078 is a selective androgen receptor modulator (SARM).¹ It induces AR-mediated transcriptional activation in PC3(AR)₂ cells ($EC_{50} = 3.99$ nM), an effect that can be inhibited by the non-steroidal AR antagonist bicalutamide (Item No. 14250).

Reference

1. Zierau, O., Kolodziejczyk, A., Vollmer, G., *et al.* Comparison of the three SARMs RAD-140, GLPG0492 and GSK-2881078 in two different *in vitro* bioassays, and in an *in silico* androgen receptor binding assay. *J. Steroid Biochem. Mol. Biol.* **189**, 81-86 (2019).

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, 11/10/2022

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