

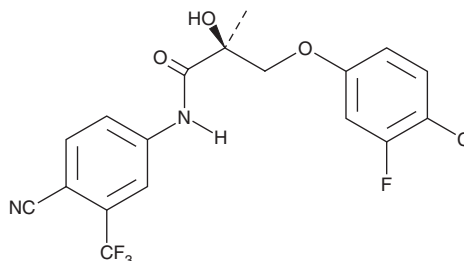
PRODUCT INFORMATION



S-23

Item No. 27304

CAS Registry No.: 1010396-29-8
Formal Name: (2S)-3-(4-chloro-3-fluorophenoxy)-N-[4-cyano-3-(trifluoromethyl)phenyl]-2-hydroxy-2-methyl-propanamide
MF: C₁₈H₁₃ClF₄N₂O₃
FW: 416.8
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 271 nm
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

S-23 is supplied as a solid. A stock solution may be made by dissolving the S-23 in the solvent of choice, which should be purged with an inert gas. S-23 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of S-23 in these solvents is approximately 10, 14, and 20 mg/ml, respectively.

S-23 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, S-23 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. S-23 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

S-23 is a selective androgen receptor modulator (SARM).¹ It binds to the AR (K_i = 1.7 nM) and induces AR-mediated transcriptional activation in CV-1 cells expressing the human receptor when used at a concentration of 10 nM. S-23 increases prostate, seminal vesicle, and levator ani muscle weights in castrated rats. It decreases testicular sperm concentration without reducing mounting behavior or the number of intromissions in intact rats when administered in combination with estradiol benzoate (Item No. 10006487). S-23 (3 mg/kg) also increases preference for sexually active intact males when administered to ovariectomized female rats.²

References

1. Jones, A., Chen, J., Hwang, D.J., *et al.* Preclinical characterization of a (S)-N-(4-cyano-3-trifluoromethylphenyl)-3-(3-fluoro, 4-chlorophenoxy)-2-hydroxy-2-methyl-propanamide: A selective androgen receptor modulator for hormonal male contraception. *Endocrinology* **150**(1), 385-395 (2009).
2. Jones, A., Hwang, D.J., Duke, C.B., III, *et al.* Nonsteroidal selective androgen receptor modulators enhance female sexual motivation. *J. Pharmacol. Exp. Ther.* **334**(2), 439-448 (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, 10/17/2022

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