

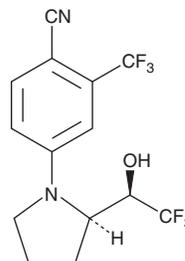
PRODUCT INFORMATION



LGD 4033

Item No. 9002046

CAS Registry No.: 1165910-22-4
Formal Name: 4-[(2R)-2-[(1R)-2,2,2-trifluoro-1-hydroxyethyl]-1-pyrrolidinyl]-2-(trifluoromethyl)-benzonitrile
MF: C₁₄H₁₂F₆N₂O
FW: 338.3
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 295 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.

Description

LGD 4033 is an orally bioavailable, nonsteroidal selective androgen receptor modulator (SARM) that binds to the androgen receptor with a K_i value of 1 nM.¹ In animal models, LGD 4033 displays robust selectivity for muscle versus prostate tissues, anabolic activity in the muscle, and anti-resorptive and anabolic activity in bone.¹ LGD 4033 has been abused as a performance-enhancing drug and the identification of its *in vitro* metabolites has been described.^{2,3} This product is intended for research and forensic applications.

References

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).
2. Geldof, L., Pozo, O.J., Lootens, L., *et al.* *In vitro* metabolism study of a black market product containing SARM LGD-4033. *Drug Test Anal.* **9(2)**, 168-178(2016).
3. Thevis, M., Lagojda, A., Kuehne, D., *et al.* Characterization of a non-approved selective androgen receptor modulator drug candidate sold via the internet and identification of *in vitro* generated phase-I metabolites for human sports drug testing. *Rapid Commun Mass Spectrom.* **29(11)**, 991-999 (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, 11/22/2022

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