

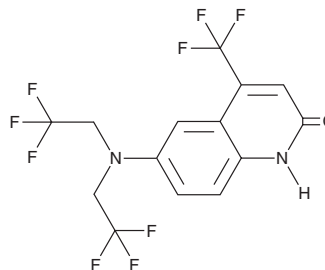
PRODUCT INFORMATION



LGD 2226

Item No. 38764

CAS Registry No.: 328947-93-9
Formal Name: 6-[bis(2,2,2-trifluoroethyl)amino]-4-(trifluoromethyl)-2(1H)-quinolinone
MF: C₁₄H₉F₉N₂O
FW: 392.2
Purity: ≥98%
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

LGD 2226 is supplied as a solid. A stock solution may be made by dissolving the LGD 2226 in the solvent of choice, which should be purged with an inert gas. LGD 2226 is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Description

LGD 2226 is an androgen receptor agonist.¹ It induces reporter activation in a reporter assay using CV-1 cells expressing human androgen receptors (EC₅₀ = 0.2 nM). LGD 2226 (10 mg/kg per day) increases prostate and levator ani muscle weight in orchidectomized (ORDX) rats.² It increases bone mineral density and breaking strength and decreases periosteal mineral apposition and bone formation rates in ORDX rats when administered at doses of 3 or 10 mg/kg per day. LGD 2226 increases mounting behavior and the number of intromissions and ejaculations in ORDX rats.

References

1. van Oeveren, A., Motamedi, M., Mani, N.S., *et al.* Discovery of 6-N,N-bis(2,2,2-trifluoroethyl)amino-4-trifluoromethylquinolin-2(1H)-one as a novel selective androgen receptor modulator. *J. Med. Chem.* **49**(21), 6143-6146 (2006).
2. Miner, J.N., Chang, W., Chapman, M.S., *et al.* An orally active selective androgen receptor modulator is efficacious on bone, muscle, and sex function with reduced impact on prostate. *Endocrinology* **148**(1), 363-373 (2007).

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, 02/13/2025

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