

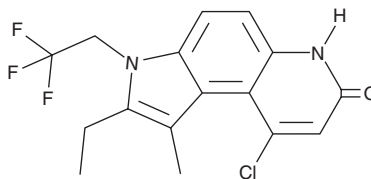
PRODUCT INFORMATION



LGD 3033

Item No. 40143

CAS Registry No.: 917891-35-1
Formal Name: 9-chloro-2-ethyl-3,6-dihydro-1-methyl-3-(2,2,2-trifluoroethyl)-7H-pyrrolo[3,2-f]quinolin-7-one
MF: C₁₆H₁₄ClF₃N₂O
FW: 342.7
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 3033 is supplied as a solid. A stock solution may be made by dissolving the LGD 3033 in the solvent of choice, which should be purged with an inert gas. LGD 3033 is soluble in acetonitrile and DMSO.

Description

LGD 3033 is an androgen receptor agonist.¹ It selectively binds to the androgen receptor ($K_i = 0.9$ nM for the human receptor) over mineralocorticoid, glucocorticoid, and progesterone receptors (K_i s = 1,261, 581, and 136 nM, respectively, for the human receptors). It also selectively induces reporter transcription in a reporter assay using CV-1 cells expressing the androgen receptor ($EC_{50} = 3.6$ nM) over CV-1 cells expressing the mineralocorticoid or progesterone receptors (EC_{50} s = 3,695 and 2,233 nM, respectively). LGD 3033 (3 mg/kg per day) increases levator ani muscle mass, but not prostate muscle mass, in orchietomized (ORDX) male rats. It increases body weight and gastrocnemius mass, bone mineral content and density in the lumbar spine and mid-femur, and bone formation rate in ovariectomized (OVX) female rats when administered at a dose of 3 mg/kg per day. LGD 3033 (30 mg/kg per day) increases the time spent in the male area in a sexual preference test by sexually experienced, but not sexually inexperienced, estradiol benzoate- and progesterone-primed OVX female rats.²

References

1. Vajda, E.G., Hogue, A., Griffiths, K.N., *et al.* Combination treatment with a selective androgen receptor modulator (SARM) and a bisphosphonate has additive effects in osteopenic female rats. *J. Bone Miner. Res.* **24**(2), 231-240 (2009).
2. Kudwa, A.E., López, F.J., McGivern, R.F., *et al.* A selective androgen receptor modulator enhances male-directed sexual preference, proceptive behavior, and lordosis behavior in sexually experienced, but not sexually naive, female rats. *Endocrinology* **151**(6), 2659-2668 (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, 02/08/2024

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