

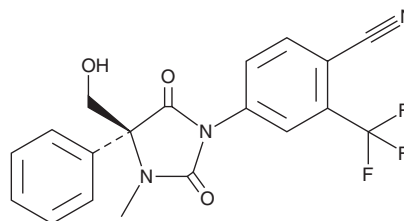
PRODUCT INFORMATION



GLPG0492

Item No. 29938

CAS Registry No.: 1215085-92-9
Formal Name: 4-[(4S)-4-(hydroxymethyl)-3-methyl-2,5-dioxo-4-phenyl-1-imidazolidinyl]-2-(trifluoromethyl)-benzonitrile
MF: C₁₉H₁₄F₃N₃O₃
FW: 389.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

GLPG0492 is supplied as a solid. A stock solution may be made by dissolving the GLPG0492 in the solvent of choice, which should be purged with an inert gas. GLPG0492 is slightly soluble in ethanol and DMSO.

Description

GLPG0492 is a selective androgen receptor modulator (SARM).¹ It activates the androgen receptor in a transactivation assay (EC₅₀ = 12 nM in HeLa cells expressing the human receptor). GLPG0492 (10 mg/kg per day) increases levator ani muscle weight, a marker of anabolic activity, but has no effect on ventral prostate weight, a marker of androgenic activity, in castrated rats. It increases twitch tension in isolated diaphragm strips, as well as increases the total distance run in a treadmill test and decreases gastrocnemius muscle fibrosis in the exercised-*mdx* mouse model of muscular dystrophy when administered at a dose of 30 mg/kg.² GLPG0492 also reverses immobilization-induced muscle atrophy in a mouse model of hindlimb immobilization.³

References

1. Nique, F., Hebbe, S., Triballeau, N., *et al.* Identification of a 4-(hydroxymethyl)diarylhydantoin as a selective androgen receptor modulator. *J. Med. Chem.* **55**(19), 8236-8247 (2012).
2. Cozzoli, A., Capogrosso, R.F., Sblendorio, V.T., *et al.* GLPG0492, a novel selective androgen receptor modulator, improves muscle performance in the exercised-*mdx* mouse model of muscular dystrophy. *Pharmacol. Res.* **72**, 9-24 (2013).
3. Blanqué, R., Lapescheux, L., Auberval, M., *et al.* Characterization of GLPG0492, a selective androgen receptor modulator, in a mouse model of hindlimb immobilization. *BMC Musculoskelet. Disord.* **15**:291 (2014).

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, 03/24/2020

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