

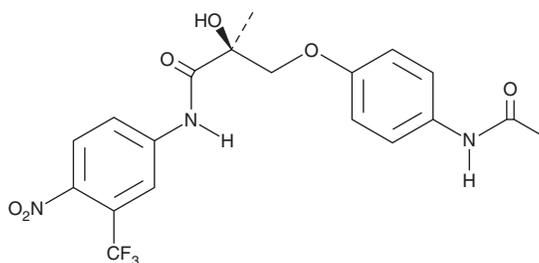
# PRODUCT INFORMATION



GTx-007

Item No. 18772

**CAS Registry No.:** 401900-40-1  
**Formal Name:** (2S)-3-[4-(acetylamino)phenoxy]-2-hydroxy-2-methyl-N-[4-nitro-3-(trifluoromethyl)phenyl]-propanamide  
**Synonyms:** Andarine, S-4, SARM S-4  
**MF:** C<sub>19</sub>H<sub>18</sub>F<sub>3</sub>N<sub>3</sub>O<sub>6</sub>  
**FW:** 441.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 248 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.

## Laboratory Procedures

GTx-007 is supplied as a crystalline solid. A stock solution may be made by dissolving the GTx-007 in the solvent of choice, which should be purged with an inert gas. GTx-007 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GTx-007 in ethanol is approximately 14 mg/ml and approximately 20 mg/ml in DMSO and DMF.

GTx-007 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GTx-007 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GTx-007 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

GTx-007 is a non-steroidal selective androgen receptor modulator (SARM) that has tissue-selective androgenic and anabolic effects in animals.<sup>1</sup> It improves muscle strength and prevents bone loss in orchidectomized rats.<sup>2</sup> Like other SARMs, GTx-007 produces anabolic effects in muscle and bone without androgenic effects.<sup>3</sup> Anabolic agents, including GTx-007 and other SARMs, are prohibited by the World Anti-Doping Agency.<sup>4</sup>

## References

1. Kearbey, J.D., Wu, D., Gao, W., *et al.* Pharmacokinetics of S-3-(4-acetylamino-phenoxy)-2-hydroxy-2-methyl-N-(4-nitro-3-trifluoromethyl-phenyl)-propionamide in rats, a non-steroidal selective androgen receptor modulator. *Xenobiotica* **34(3)**, 273-280 (2004).
2. Gao, W., Reiser, P.J., Coss, C.C., *et al.* Selective androgen receptor modulator treatment improves muscle strength and body composition and prevents bone loss in orchidectomized rats. *Endocrinology* **146(11)**, 4887-4897 (2005).
3. Girgis, C.M., Mokbel, N., and DiGirolamo, D.J. Therapies for musculoskeletal disease: Can we treat two birds with one stone? *Curr. Osteoporos. Rep.* **12(2)**, 142-153 (2014).
4. Geyer, H., Schänzner, W., and Thevis, M. Anabolic agents: Recent strategies for their detection and protection from inadvertent doping. *Br. J. Sports Med.* **48(10)**, 820-826 (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, 11/28/2022

## 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