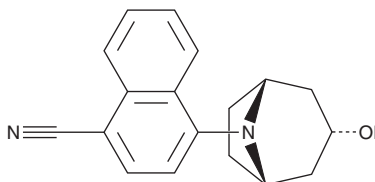


# PRODUCT INFORMATION



**AC-262536**  
Item No. 31598

**CAS Registry No.:** 870888-46-3  
**Formal Name:** 4-[(3-*endo*)-3-hydroxy-8-azabicyclo[3.2.1]oct-8-yl]-1-naphthalenecarbonitrile  
**MF:** C<sub>18</sub>H<sub>18</sub>N<sub>2</sub>O  
**FW:** 278.4  
**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

AC-262536 is supplied as a solid. A stock solution may be made by dissolving the AC-262536 in the solvent of choice, which should be purged with an inert gas. AC-262536 is soluble in DMSO.

## Description

AC-262536 is a selective androgen receptor modulator (SARM; K<sub>i</sub> = 5.01 nM).<sup>1</sup> It is also a partial agonist of the androgen receptor (EC<sub>50</sub> = 1.58 nM in a luciferase assay). AC-262536 is selective for the androgen receptor over a panel of 47 additional human nuclear receptors at 10 μM. It inhibits dihydroxytestosterone-induced proliferation of LNCaP prostate cancer cells when used at concentrations of 0.1 and 1 μM. AC-262536 (3, 10, and 30 mg/kg) reduces plasma luteinizing hormone levels and increases levator ani muscle weight, a marker of anabolic activity, in castrated rats. It exhibits smaller increases in prostate gland and seminal vesicle weight, markers of androgenic activity, compared with testosterone in the same model.

## Reference

1. Piu, F., Gardell, L.R., Son, T., *et al.* Pharmacological characterization of AC-262536, a novel selective androgen receptor modulator. *J. Steroid Biochem. Mol. Biol.* **109**(1-2), 129-137 (2008).

**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, 09/11/2020

## 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