

PRODUCT INFORMATION



Ibutamoren

Item No. 44475

CAS Registry No.: 159634-47-6
Formal Name: 2-amino-N-[(1R)-2-[1,2-dihydro-1-(methylsulfonyl)spiro[3H-indole-3,4'-piperidin]-1'-yl]-2-oxo-1-[(phenylmethoxy)methyl]ethyl]-2-methyl-propanamide

Synonyms: L-163,191, MK-0677

MF: C₂₇H₃₆N₄O₅S

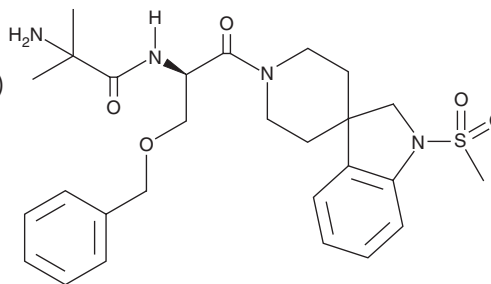
FW: 528.7

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.

Description

Ibutamoren (Item No. 44475) is an analytical reference standard categorized as a growth hormone secretagogue.¹ Ibutamoren has been detected in urine from competition athletes.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 18003).

References

1. Patchett, A.A., Nargund, R.P., Tata, J.R., *et al.* Design and biological activities of L-163,191 (MK-0677): A potent, orally active growth hormone secretagogue. *Proc. Natl. Acad. Sci. USA* **92**(15), 7001-7005 (1995).
2. Börjesson, A., Lehtihet, M., Andersson, A., *et al.* Studies of athlete biological passport biomarkers and clinical parameters in male and female users of anabolic androgenic steroids and other doping agents. *Drug Test Anal.* **12**(4), 514-523 (2020).

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, 10/03/2025

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