

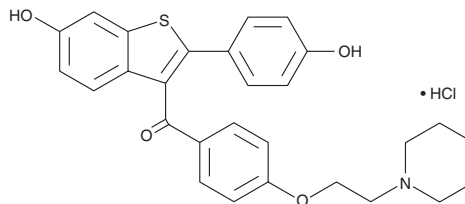
PRODUCT INFORMATION



Raloxifene (hydrochloride)

Item No. 27189

CAS Registry No.: 82640-04-8
Formal Name: [6-hydroxy-2-(4-hydroxyphenyl)benzo[b]thien-3-yl][4-[2-(1-piperidinyl)ethoxy]phenyl]-methanone, monohydrochloride
Synonyms: Keoxifene, LY156758
MF: C₂₈H₂₇NO₄S • HCl
FW: 510.0
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 282 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Raloxifene (hydrochloride) (Item No. 27189) is an analytical reference standard categorized as a selective estrogen receptor modulator (SERM).¹ It has been detected as an adverse analytical finding (AAF) during anti-doping testing.² This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 10011620).

References

1. Sun, D., Jones, N.R., Manni, A., *et al.* Characterization of raloxifene glucuronidation: Potential role of UGT1A8 genotype on raloxifene metabolism *in vivo*. *Cancer Prev. Res. (Phila)*. **6(7)**, 719-730 (2013).
2. World Anti-Doping Agency. *2014 Anti-doping testing figures report* (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, 07/25/2023

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