

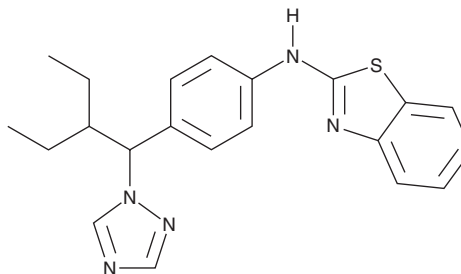
PRODUCT INFORMATION



Talarozole

Item No. 42372

CAS Registry No.: 201410-53-9
Formal Name: N-[4-[2-ethyl-1-(1H-1,2,4-triazol-1-yl)butyl]phenyl]-2-benzothiazolamine
Synonyms: R115866, Rambazole
MF: C₂₁H₂₃N₅S
FW: 377.5
Purity: ≥95%
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.

Laboratory Procedures

Talarozole is supplied as a solid. A stock solution may be made by dissolving the talarozole in the solvent of choice, which should be purged with an inert gas. Talarozole is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

Talarozole is an inhibitor of cytochrome P450 (CYP) isoform 26 (CYP26; IC₅₀ = 0.004 μM).¹ It is selective for CYP26 over CYP2A1, CYP2C11, CYP17, and CYP19 (IC₅₀s = 9.1, 7.2, 2.6, and 1.3 μM, respectively). Talarozole (2.5 mg/kg) increases the levels of all-*trans* retinoic acid in plasma, skin, liver, fat, kidneys, and spleen in rats. It inhibits estrogen-induced vaginal keratinization in ovariectomized rats (ED₅₀ = 1 mg/kg), an effect that can be reversed by AGN 193109 (Item No. 23975). Talarozole (0.6-2.6 mg/kg) increases pinnal epidermal thickness and tail epidermal orthokeratosis percentage in mice.

Reference

1. Stoppie, P., Borgers, M., Borghgraef, P., *et al.* R115866 inhibits all-*trans*-retinoic acid metabolism and exerts retinoidal effects in rodents. *J. Pharmacol. Exp. Ther.* **293**(1), 304-312 (2000).

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, 12/16/2024

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