

PRODUCT INFORMATION



ZK 164015

Item No. 45945

CAS Registry No.: 177583-70-9

Formal Name: 2-(4-hydroxyphenyl)-3-methyl-1-[10-(pentylsulfonyl)decyl]-1H-indol-5-ol

MF: C₃₀H₄₃NO₄S

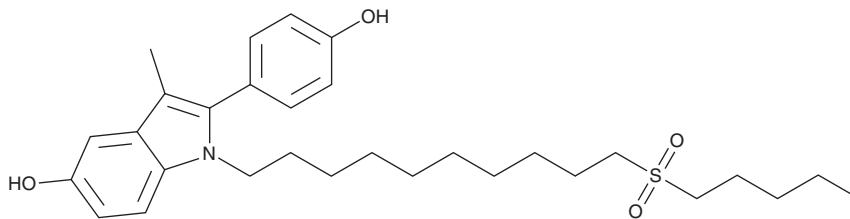
FW: 513.7

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

ZK 164015 is supplied as a solid. A stock solution may be made by dissolving the ZK 164015 in the solvent of choice, which should be purged with an inert gas. ZK 164015 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

ZK 164015 is an estrogen receptor antagonist (IC₅₀ = 25 nM in a reporter assay).¹ It inhibits the proliferation of MCF-7 breast cancer cells stimulated with E₂ (17β-estradiol; IC₅₀ = 51 nM) and unstimulated MCF-7 cells (IC₅₀ = 1.4 nM). ZK 164015 does not increase uterine weight in mice.² It also inhibits the growth of *P. falciparum* in isolated human red blood cells in a concentration-dependent manner.³

References

1. Walter, G., Liebl, R. and von Angerer, E. Synthesis and biological evaluation of stilbene-based pure estrogen antagonists. *Bioorg. Med. Chem. Lett.* **14**(18), 4659-4663 (2004).
2. Biberger, C. and von Angerer, E. 2-Phenylindoles with sulfur containing side chains. Estrogen receptor affinity, antiestrogenic potency, and antitumor activity. *J. Steroid Biochem. Mol. Biol.* **58**(1), 31-43 (1996).
3. Honfozo, A., Hougue, R., Vandeputte, A., et al. An image-based high-content screening for compounds targeting *Toxoplasma gondii* repurposed inhibitors effective against the malaria parasite *Plasmodium falciparum*. *Front. Cell. Infect. Microbiol.* **13**, 1102551 (2023).

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, 07/01/2026

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