

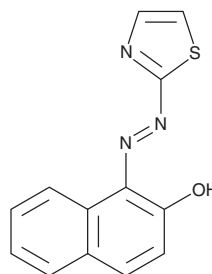
PRODUCT INFORMATION



ERGi-USU

Item No. 30875

CAS Registry No.: 1147-56-4
Formal Name: 1-[2-(2-thiazolyl)diazenyl]-2-naphthalenol
Synonym: NSC 139021
MF: C₁₃H₉N₃OS
FW: 255.3
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 226, 482 nm
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

ERGi-USU is supplied as a solid. A stock solution may be made by dissolving the ERGi-USU in the solvent of choice, which should be purged with an inert gas. ERGi-USU is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ERGi-USU in these solvents is approximately 0.5, 5, and 15 mg/ml, respectively.

Description

ERGi-USU is an inhibitor of the oncoprotein ETS-related gene (*ERG*; IC₅₀ = 315 nM).¹ It inhibits proliferation of *ERG*-expressing VCaP prostate, COLO 320 colon, and KG-1 and MOLT-4 leukemia cancer cells (IC₅₀s = 30-400 nM) but not LNCaP, LAPC4, and MDA PCa 2b prostate cancer and human umbilical vein endothelial cells (HUVECs), which do not express *ERG* (IC₅₀s = >10 μM for all). ERGi-USU (100 and 150 mg/kg) reduces tumor growth in a VCaP mouse xenograft model.

Reference

1. Mohamed, A.A., Xavier, C.P., Sukumar, G., *et al.* Identification of a small molecule that selectively inhibits ERG-positive cancer cell growth. *Cancer Res.* **78(13)**, 3659-3671 (2018).

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, 11/30/2022

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