

PRODUCT INFORMATION

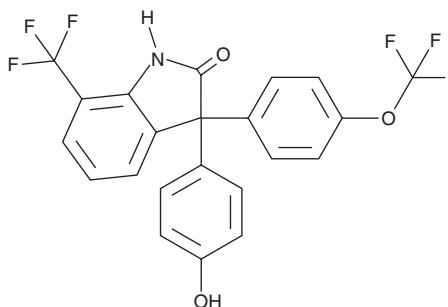


(±)-ErSO

Item No. 39676

CAS Registry No.: 2407860-40-4
Formal Name: N-[[[(1S)-3-carboxy-1-[(1,1-dimethylethoxy)carbonyl]propyl]amino]carbonyl]-L-glutamic acid, 1,5-bis(1,1-dimethylethyl) ester

MF: C₂₂H₁₃F₆NO₃
FW: 453.3
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

(±)-ErSO is supplied as a solid. A stock solution may be made by dissolving the (±)-ErSO in the solvent of choice, which should be purged with an inert gas. (±)-ErSO is soluble in dimethyl formamide and acetonitrile. (±)-ErSO is slightly soluble in DMSO and ethanol.

Description

(±)-ErSO is an activator of the anticipatory unfolded protein response (a-UPR) and a racemic mixture of the a-UPR activator (R)-ErSO (Item No. 35684) and its inactive enantiomer (S)-ErSO.¹ (±)-ErSO is cytotoxic to MCF-7 cells (IC₅₀ = 42.6 nM), as well as T47D cells when used at concentrations of 0.5 and 1 μM.

Reference

1. Boudreau, M.W., Duraki, D., Wang, L., *et al.* A small-molecule activator of the unfolded protein response eradicates human breast tumors in mice. *Sci. Transl. Med.* **13(603)**, eabf1383 (2021).

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, 02/08/2024

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