

PRODUCT INFORMATION

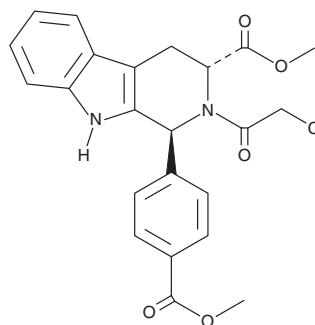


(1S,3R)-RSL3

Item No. 19288

CAS Registry No.: 1219810-16-8
Formal Name: (1S,3R)-2-(2-chloroacetyl)-2,3,4,9-tetrahydro-1-[4-(methoxycarbonyl)phenyl]-1H-pyrido[3,4-b]indole-3-carboxylic acid, methyl ester

MF: C₂₃H₂₁ClN₂O₅
FW: 440.9
Purity: ≥98%
UV/Vis.: λ_{max}: 220 nm
Supplied as: A crystalline 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

(1S,3R)-RSL3 is supplied as a crystalline solid. A stock solution may be made by dissolving the (1S,3R)-RSL3 in the solvent of choice, which should be purged with an inert gas. (1S,3R)-RSL3 is soluble in organic solvents such as ethanol, DMSO and dimethyl formamide (DMF). The solubility of (1S,3R)-RSL3 in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

(1S,3R)-RSL3 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (1S,3R)-RSL3 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (1S,3R)-RSL3 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(1S,3R)-RSL3 is a ferroptosis inducer and is canonically known as an inhibitor of glutathione peroxidase 4 (GPX4).¹⁻³ It inhibits GPX4 in a context-dependent manner: (1S,3R)-RSL3 inhibits GPX4 in the presence of cell cytosol or recombinant human 14-3-3ε, an adaptor protein, but does not inhibit recombinant human selenocysteine-containing GPX4 or purified rat testis GPX4, when used at a concentration of 100 μM.⁴⁻⁵ (1S,3R)-RSL3 (50 nM) inhibits peroxidase activity in GPX4-overexpressing COH-BR1 breast cancer cell lysates.¹ It also inhibits thioredoxin reductase 1 (TrxR1; IC₅₀ = 7.9 μM in a cell-free assay).⁴ (1S,3R)-RSL3 is selectively cytotoxic to H-Ras^{V12}-expressing BJeLR cells over wild-type Ras-expressing BJeH cells (EC₅₀s = 0.01 and 2 μM, respectively) and induces lipid peroxidation, a hallmark of ferroptosis, in Pfa1 fibroblasts when used at a concentration of 100 nM.^{1,2} *In vivo*, (1S,3R)-RSL3 (100 mg/kg) prevents tumor formation and inhibits tumor growth in BJeLR mouse xenograft models.¹

References

1. Yang, W.S., SriRamaratnam, R., Welsch, M.E., *et al.* Regulation of ferroptotic cancer cell death by GPX4. *Cell* **156**(1-2), 317-331 (2015).
2. Kagan, V.E., Mao, G., Qu, F., *et al.* Oxidized arachidonic and adrenic PEs navigate cells to ferroptosis. *Nat. Chem. Biol.* **13**(1), 81-90 (2017).
3. Yang, W.S. and Stockwell, B.R. Ferroptosis: Death by lipid peroxidation. *Trends Cell Biol.* **26**(3), 165-176 (2016).
4. Cheff, D.M., Huang, C., Scholzen, K.C., *et al.* The ferroptosis inducing compounds RSL3 and ML162 are not direct inhibitors of GPX4 but of TXNRD1. *Redox Biol.* **62**:102703, (2023).
5. Vučković, A.-M., Travain, V.B., Bordin, L., *et al.* Inactivation of the glutathione peroxidase GPx4 by the ferroptosis-inducing molecule RSL3 requires the adaptor protein 14-3-3ε. *FEBS Lett.* **594**(4), 611-624 (2020).

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, 06/21/2023

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