

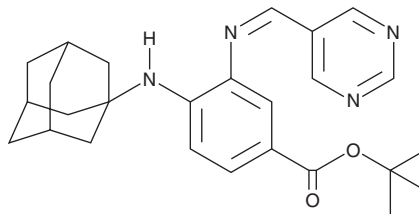
PRODUCT INFORMATION



SRS16-86

Item No. 26752

CAS Registry No.: 1793052-96-6
Formal Name: 3-[(Z)-(5-pyrimidinylmethylene)amino]-4-(tricyclo[3.3.1.1^{3,7}]dec-1-ylamino)-benzoic acid, 1,1-dimethylethyl ester
MF: C₂₆H₃₂N₄O₂
FW: 432.6
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 303, 429 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

SRS16-86 is supplied as a crystalline solid. A stock solution may be made by dissolving the SRS16-86 in the solvent of choice. SRS16-86 is soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 25 mg/ml. SRS16-86 is also slightly soluble in ethanol and DMSO.

Description

SRS16-86 is an inhibitor of ferroptosis.¹ It inhibits ferroptosis induced by erastin (Item No. 17754) in HT-1080 and NIH3T3 cells when used at a concentration of 1 μM. SRS16-86 (2 mg/kg) prevents renal tubular damage and increases in serum levels of urea and creatine in a mouse model of renal ischemia-reperfusion injury (IRI). In a rat model of spinal cord injury, SRS16-86 (15 mg/kg) increases the levels of glutathione peroxidase 4 (GPX4), system x_c⁻ cystine/glutamate transporter (xCT), and glutathione (GSH) and decreases levels of IL-1β, TNF-α, ICAM-1, and the lipid peroxidation marker 4-hydroxy nonenal (4HNE) in injured spinal cord epicenters.² It also increases tissue sparing and improves locomotor recovery in the same model.

References

- Linkermann, A., Skouta, R., Himmerkus, N., *et al.* Synchronized renal tubular cell death involves ferroptosis. *Proc. Natl. Acad. Sci. U.S.A.* **111**(47), 16836-16841 (2014).
- Zhang, Y., Sun, C., Zhao, C., *et al.* Ferroptosis inhibitor SRS 16-86 attenuates ferroptosis and promotes functional recovery in contusion spinal cord injury. *Brain Res.* **1706**, 48-57 (2019).

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, 02/16/2024

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