

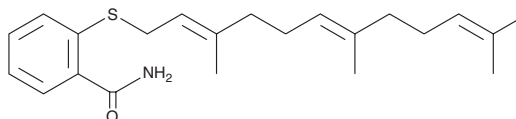
PRODUCT INFORMATION



Farnesyl Thiosalicylic Acid Amide

Item No. 13175

CAS Registry No.: 1092521-74-8
Formal Name: 2-[[[(2E,6E)-3,7,11-trimethyl-2,6,10-dodecatrien-1-yl]thio]-benzamide
Synonyms: FTS Amide, Salirasib Amide
MF: C₂₂H₃₁NOS
FW: 357.6
Purity: ≥96%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Farnesyl thiosalicylic acid (FTS) amide is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of FTS amide in these solvents is approximately 10 mg/ml.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Association of Ras protein with the inner surface of the plasma membrane is required for Ras signaling activity. FTS is an inhibitor of Ras-mediated signaling that functions by dislodging Ras from the cell membrane thereby rendering it susceptible to proteolytic degradation.¹ FTS amide is an FTS derivative with an amide added to the carboxyl group. FTS amide inhibits the growth of PANC-1 and U87 tumor cells with IC₅₀ values of 20 and 10 μM, respectively, a relatively higher potency compared to that of FTS (IC₅₀s = 35 and 50 μM, respectively).² Treatment of nude mice bearing either U87 glioblastoma or PANC-1 tumors with 100 mg/kg FTS-amide twice daily for four days inhibits tumor growth by at least 50% of controls.

References

1. Haklai, R., Weisz, M.G., Elad, G., *et al.* Dislodgment and accelerated degradation of ras. *Biochemistry* **37**, 1306-1314 (1998).
2. Goldberg, L., Haklai, R., Bauer, V., *et al.* New derivatives of farnesylthiosalicylic acid (salirasib) for cancer treatment: Farnesylthiosalicylamide inhibits tumor growth in nude mice models. *J. Med. Chem.* **52**, 197-205 (2009).

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, 04/05/2024

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