

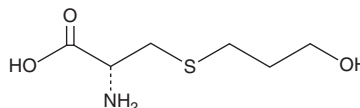
PRODUCT INFORMATION



Fudosteine

Item No. 16722

CAS Registry No.: 13189-98-5
Formal Name: S-(3-hydroxypropyl)-L-cysteine
Synonym: SS320A
MF: C₆H₁₃NO₃S
FW: 179.2
Purity: ≥95%
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

Fudosteine is supplied as a crystalline solid. Aqueous solutions of fudosteine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of fudosteine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Fudosteine is a cysteine derivative that interferes with the increase in goblet cell number, expression of the mucin MUC5AC, and subsequent mucin secretion, in airways agonized with tobacco smoke, isoproterenol, lipopolysaccharide, TNF- α , or oxidants.¹⁻³ Fudosteine, which is a thiol compound with antioxidant properties, limits NF- κ B signaling and reduces inflammatory gene expression.³ It has greater bioavailability than N-acetyl-cysteine and increases cysteine levels in cells.³

References

1. Takahashi, K., Mizuno, H., Ohno, H., *et al.* Effects of SS320A, a new cysteine derivative, on the change in the number of goblet cells induced by isoproterenol in rat tracheal epithelium. *Jpn. J. Pharmacol.* **77(1)**, 71-77 (1998).
2. Rhee, C.K., Kang, C.M., You, M.B., *et al.* Effect of fudosteine on mucin production. *Eur. Respir. J.* **32(5)**, 1195-1202 (2008).
3. Rahman, I. Pharmacological antioxidant strategies as therapeutic interventions for COPD. *Biochim. Biophys. Acta* **1822(5)**, 714-728 (2012).

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, 11/09/2022

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