

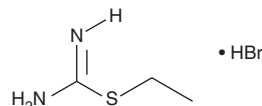
PRODUCT INFORMATION



S-ethyl Isothiourea (hydrobromide)

Item No. 81275

CAS Registry No.: 1071-37-0
Formal Name: ethyl carbamimidothioate, hydrobromide
Synonyms: Ethiron bromide; 2-Ethyl-2-thiopseudourea hydrobromide; SEIT (HBr); WR 539
MF: C₃H₈N₂S • HBr
FW: 185.1
Purity: ≥98%
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

S-ethyl Isothiourea (hydrobromide) (SEIT (HBr)) is supplied as a crystalline solid. A stock solution may be made by dissolving the SEIT (HBr) in an organic solvent purged with an inert gas. SEIT (HBr) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SEIT (HBr) in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of SEIT (HBr) can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of SEIT (HBr) in PBS, (pH 7.2), is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

SEIT is a potent inhibitor of NOS *in vitro* but lacks good *in vivo* efficacy due to poor cellular penetration. The K_i values are 19, 39 and 29 nM using purified human iNOS, eNOS, and nNOS, respectively.¹⁻³

References

1. Garvey, E.P., Oplinger, J.A., Tanoury, G.J., *et al.* Potent and selective inhibition of human nitric oxide synthases. Inhibition by non-amino acid isothioureas. *J. Biol. Chem.* **269**, 26669-26676 (1994).
2. Nakane, M., Klinghofer, V., Kuk, J.E., *et al.* Novel potent and selective inhibitors of inducible nitric oxide synthase. *Mol. Pharmacol.* **47**, 831-834 (1995).
3. Southan, G.J., Szabó, C., and Thiemermann, C. Isothioureas: Potent inhibitors of nitric oxide synthases with variable isoform selectivity. *Br. J. Pharmacol.* **114**, 510-516 (1995).

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, 11/17/2022

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