

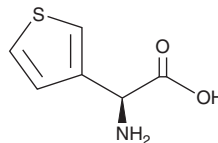
PRODUCT INFORMATION



(S)-3-Thienylglycine

Item No. 30912

CAS Registry No.: 1194-87-2
Formal Name: αS-amino-3-thiopheneacetic acid
Synonyms: L-R-(3-Thienyl)glycine, L-α-3-Thienylglycine
MF: C₆H₇NO₂S
FW: 157.2
Purity: ≥95%
UV/Vis.: λ_{max}: 233 nm
Supplied as: A 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)-3-Thienylglycine is supplied as a solid. A stock solution may be made by dissolving the (S)-3-thienylglycine in water. (S)-3-Thienylglycine is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

(S)-3-Thienylglycine is a thienyl-containing amino acid.¹ It has been used as an affinity ligand to adsorb IgM and IgG antibodies from isolated human plasma.

Reference

1. Hatanaka, Y., Tsukiji, M., Itoh, A., *et al.* Selective adsorption of immunoglobulin G and immunoglobulin M from plasma without adsorption of fibrinogen by using thienyl amino acids as ligands. *J. Chromatogr.* **4(6)**, 1000190 (2013).

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, 11/04/2022

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