

PRODUCT INFORMATION



Somatostatin-28 (1-14) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt)

Item No. 24992

Formal Name: L-seryl-L-alanyl-L-asparaginyl-L-seryl-L-asparaginyl-L-prolyl-L-alanyl-L-methionyl-L-alanyl-L-prolyl-L-arginyl-L- α -glutamyl-L-arginyl-L-lysine, trifluoroacetate salt
MF: $C_{61}H_{105}N_{23}O_{21}S \cdot XCF_3COOH$
FW: 1,528.7
Purity: $\geq 95\%$
Supplied as: A lyophilized powder
Storage: $-20^\circ C$
Stability: ≥ 4 years

H—Ser—Ala—Asn—Ser—Asn—Pro—Ala—Met—Ala—Pro—Arg—Glu—Arg—Lys—OH
• XCF₃COOH

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

Laboratory Procedures

Somatostatin-28 (1-14) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the somatostatin-28 (1-14) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) in water. The solubility of somatostatin-28 (1-14) (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Somatostatin-28 (1-14) is an N-terminal fragment of the neuropeptide somatostatin-28 (Item No. 24991).¹ It has been used to produce antibodies selective for somatostatin-28 that do not bind to the C-terminal fragment somatostatin-14 (Item No. 20809).²

References

1. He, H.T., Johnson, K., Thermos, K., *et al.* Purification of a putative brain somatostatin receptor. *Proc. Natl. Acad. Sci. USA* **86(5)**, 1480-1484 (1989).
2. Zeggari, M., Esteve, J.P., Raully, I., *et al.* Co-purification of a protein tyrosine phosphatase with activated somatostatin receptors from rat pancreatic acinar membranes. *Biochem. J.* **303(Pt 2)**, 441-448 (1994).

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, 11/18/2022

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