

PRODUCT INFORMATION



Crosstide (trifluoroacetate salt)

Item No. 36889

Formal Name: glycyl-L-arginyl-L-prolyl-L-arginyl-L-threonyl-L-seryl-L-seryl-L-phenylalanyl-L-alanyl-L-α-glutamyl-glycine, trifluoroacetate salt
Synonyms: Gly-Arg-Pro-Arg-Thr-Ser-Ser-Phe-Ala-Glu-Gly-OH, GRPRTSSFAEG-OH

MF: C₄₈H₇₇N₁₇O₁₇ • XCF₃COOH

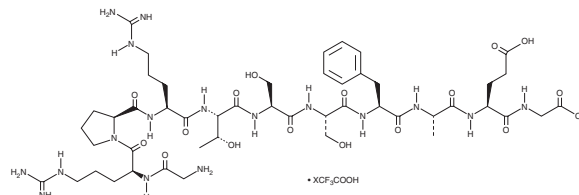
FW: 1,164.2

Purity: ≥95%

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

Crosstide (trifluoroacetate salt) is supplied as a solid. Aqueous solutions of crosstide (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of crosstide (trifluoroacetate salt) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Crosstide is a synthetic peptide substrate for Akt.¹ It corresponds to the amino acid sequence of glycogen synthase kinase 3 (GSK3) that surrounds the serine residue phosphorylated by p90 ribosomal S6 kinase 1 (RSK1) or p70 ribosomal S6 kinase (p70S6K). Crosstide has been used in the study of Akt activity.^{1,2} It is also a substrate for serum/glucocorticoid-regulated kinase family member 3 (SGK3).³

References

1. Cross, D.A.E., Slessi, D.R., Cohen, P., *et al.* Inhibition of glycogen synthase kinase-3 by insulin mediated by protein kinase B. *Nature* **378**(6559), 785-789 (1995).
2. Tanti, J.-F., Grillo, S., Gremeaux, T., *et al.* Potential role of protein kinase B in glucose transporter 4 translocation in adipocytes. *Endocrinology* **138**(5), 2005-2010 (1997).
3. Dai, F., Yu, L., He, H., *et al.* Human serum and glucocorticoid-inducible kinase-like kinase (SGKL) phosphorylates glycogen synthase kinase 3 beta (GSK-3β) at serine-9 through direct interaction. *Biochem. Biophys. Res. Commun.* **293**(4), 1191-1196 (2002).

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, 12/21/2022

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