

PRODUCT INFORMATION



Autocamtide 2 (trifluoroacetate salt)

Item No. 24788

Formal Name: L-lysyl-L-lysyl-L-alanyl-L-leucyl-L-arginyl-L-arginyl-L-glutaminy-L- α -glutamyl-L-threonyl-L-valyl-L- α -aspartyl-L-alanyl-L-leucine, trifluoroacetate salt
MF: $C_{65}H_{118}N_{22}O_{20} \cdot XCF_3COOH$
FW: 1,527.8
Purity: $\geq 95\%$
Supplied as: A lyophilized powder
Storage: $-20^\circ C$
Stability: ≥ 4 years

H-Lys-Lys-Ala-Leu-Arg-Arg-Gln-Glu-Thr-Val-Asp-Ala-Leu-OH
 $\cdot XCF_3COOH$

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

Laboratory Procedures

Autocamtide 2 (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the autocamtide 2 (trifluoroacetate salt) in water. The solubility of autocamtide 2 (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Autocamtide 2 is a peptide substrate for calcium/calmodulin-dependent protein kinase (CaMK) and CaMKII.^{1,2} It is selectively phosphorylated by CaMK over bovine brain protein kinase C (PKC; K_m s = 2 and $>50 \mu M$, respectively).² Autocamtide 2 phosphorylation can be used to quantify CaMK and CaMKII activity.

References

- Hanson, P.I., Kapiloff, M.S., Lou, L.L., *et al.* Expression of a multifunctional Ca^{2+} /calmodulin-dependent protein kinase and mutational analysis of its autoregulation. *Neuron* **3(1)**, 59-70 (1989).
- Wenham, R.M., Landt, M., and Easom, R.A. Glucose activates the multifunctional Ca^{2+} /calmodulin-dependent protein kinase II in isolated rat pancreatic islets. *J. Biol. Chem.* **269(7)**, 4947-4952 (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