

PRODUCT INFORMATION



Ziptide (trifluoroacetate salt)

Item No. 37466

Formal Name: L-lysyl-L-lysyl-L-leucyl-L-asparaginyl-L-arginyl-L-threonyl-L-leucyl-L-seryl-L-phenylalanyl-L-alanyl-L- α -glutamyl-L-prolyl-glycine, trifluoroacetate salt

Peptide Sequence: KKLNRRTLSFAEPG-OH

MF: $C_{65}H_{109}N_{19}O_{19} \cdot XCF_3COOH$

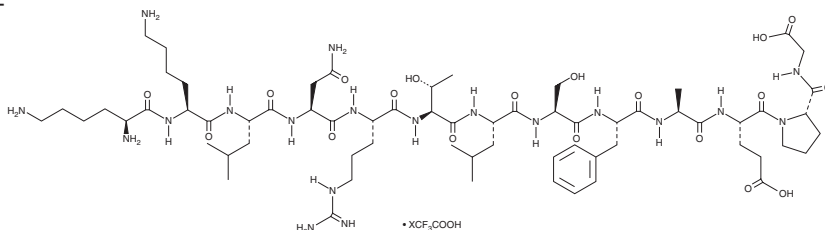
FW: 1,460.7

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Ziptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the ziptide (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Ziptide is a peptide substrate for various serine/threonine protein kinases, including MAPK activated protein kinase 2 (MAPKAPK2), MAPKAPK3, MAPKAPK5, checkpoint kinase 1 (Chk1), AMP-activated protein kinase (AMPK), and calcium/calmodulin-dependent protein kinase II (CamKII).¹ It has been used to evaluate CamKII activity for inhibitor development.²

References

1. Ross, H., Armstrong, C.G., and Cohen, P. A non-radioactive method for the assay of many serine/threonine-specific protein kinases. *Biochem J.* **366(Pt 3)**, 977-981 (2002).
2. Komiya, M., Asano, S., Koike, N., *et al.* Synthesis and structure based optimization of 2-(4-phenoxybenzoyl)-5-hydroxyindole as a novel CaMKII inhibitor. *Bioorg. Med. Chem.* **20(23)**, 6840-6847 (2012).

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, 10/05/2023

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