

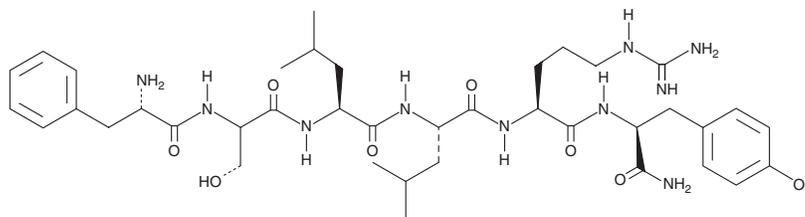
PRODUCT INFORMATION



FSLLRY-amide

Item No. 14709

CAS Registry No.: 245329-02-6
Formal Name: L-phenylalanyl-L-seryl-L-leucyl-L-leucyl-L-arginyl-L-tyrosinamide
MF: C₃₉H₆₀N₁₀O₈
FW: 797.0
Purity: ≥95%
UV/Vis.: λ_{max}: 280 nm
Supplied as: A crystalline 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

FSLLRY-amide is supplied as a crystalline solid. A stock solution may be made by dissolving the FSLLRY-amide in the solvent of choice. FSLLRY-amide is soluble in organic solvents such as acetonitrile and DMSO, which should be purged with an inert gas. The solubility of FSLLRY-amide in these solvents is approximately 0.5 and 20 mg/ml, respectively.

Description

FSLLRY-amide is a peptide antagonist of protease-activated receptor (PAR) 2 which blocks activation by trypsin (IC₅₀ = ~50 μM) but not by the synthetic agonist SLIGRL-NH₂ (Item No. 16723).¹ It does not affect thrombin-induced activation of PAR1 or trypsin-mediated proteolysis of casein.¹ It can be used to study PAR2 action in cells and in animals.^{2,3}

References

1. Al-Ani, B., Saifeddine, M., Wijesuriya, S.J., *et al.* Modified proteinase-activated receptor-1 and -2 derived peptides inhibit proteinase-activated receptor-2 activation by trypsin. *J. Pharmacol. Exp. Ther.* **300(2)**, 702-708 (2002).
2. McLarty, J.L., Meléndez, G.C., Brower, G.L., *et al.* Tryptase/protease-activated receptor 2 interactions induce selective mitogen-activated protein kinase signaling and collagen synthesis by cardiac fibroblasts. *Hypertension* **58(2)**, 264-270 (2011).
3. Chen, Y., Yang, C., and Wang, Z.J. Proteinase-activated receptor 2 sensitizes transient receptor potential vanilloid 1, transient receptor potential vanilloid 4, and transient receptor potential ankyrin 1 in paclitaxel-induced neuropathic pain. *Neuroscience* **193**, 440-451 (2011).

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, 11/23/2022

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