

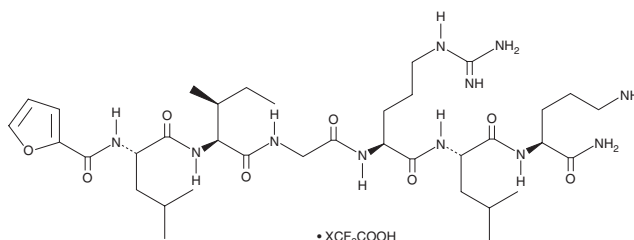
PRODUCT INFORMATION



2-furoyl-LIGRLO amide (trifluoroacetate salt)

Item No. 14594

Formal Name: N-(2-furanylcarbonyl)-L-leucyl-L-isoleucylglycyl-L-arginyl-L-leucyl-L-ornithinamide, trifluoroacetate
MF: C₃₆H₆₃N₁₁O₈ • XCF₃COOH
FW: 778.0
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

2-furoyl-LIGRLO amide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the 2-furoyl-LIGRLO amide (trifluoroacetate salt) in the solvent of choice. 2-furoyl-LIGRLO amide (trifluoroacetate salt) is soluble in water at a concentration of approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-furoyl-LIGRLO amide is a potent and selective agonist of the proteinase-activated receptor, PAR2 (pD₂ = 7.0).¹ It does not activate PAR1 or PAR4. Through its effects on PAR2, 2-furoyl-LIGRLO amide produces a concentration-dependent relaxation of arteries previously contracted with α₁ adrenergic receptor agonists.¹

Reference

- McGuire, J.J., Saifeddine, M., Triggle, C.R., *et al.* 2-Furoyl-LIGRLO-amide: A potent and selective proteinase-activated receptor 2 agonist. *J. Pharmacol. Exp. Ther.* **309**(3), 1124-1131 (2004).

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, 02/21/2019

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