

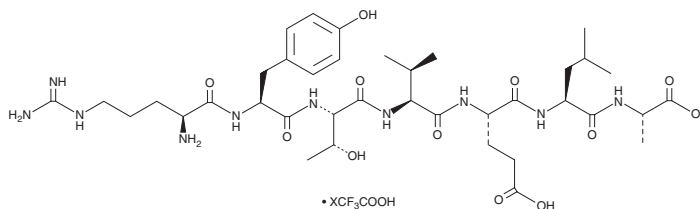
PRODUCT INFORMATION



RYTVELA (trifluoroacetate salt)

Item No. 37490

CAS Registry No.: 1200829-06-6
Formal Name: L-arginyl-L-tyrosyl-L-threonyl-L-valyl-L- α -glutamyl-L-leucyl-L-alanine
Synonym: H-Arg-Tyr-Thr-Val-Glu-Leu-Ala-OH
Peptide Sequence: RYTVELA-OH
MF: C₃₈H₆₂N₁₀O₁₂ • XCF₃COOH
FW: 851.0
Purity: ≥98%
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

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

Description

RYTVELA is a derivative of the peptide antagonist of interleukin-1 receptor 1 (IL-1R1) rytvela that contains all L-amino acids.¹

Reference

1. Quiniou, C., Kooli, E., Joyal, J.S., *et al.* Interleukin-1 and ischemic brain injury in the newborn: Development of a small molecule inhibitor of IL-1 receptor. *Semin. Perinatol.* **32(5)**, 352-333 (2008).

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, 10/30/2023

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