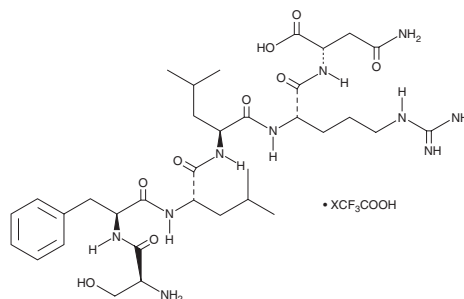


PRODUCT INFORMATION

TRAP-6 Peptide (trifluoroacetate salt)

Item No. 24125

Formal Name: L-seryl-L-phenylalanyl-L-leucyl-L-leucyl-L-arginyl-L-asparagine, trifluoroacetate salt
Synonyms: Ser-Phe-Leu-Leu-Arg-Asn, Thrombin Receptor Agonist Peptide-6
MF: $C_{34}H_{56}N_{10}O_9 \cdot XCF_3COOH$
FW: 748.9
Purity: $\geq 95\%$
Supplied as: A lyophilized powder
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

TRAP-6 peptide (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the TRAP-6 peptide (trifluoroacetate salt) in water. The solubility of TRAP-6 peptide (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

TRAP-6 peptide is a hexapeptide corresponding to residues 42-47 of protease-activated receptor 1 (PAR1).¹ It acts as an agonist of PAR1, inducing platelet aggregation in human platelet-rich plasma *ex vivo* ($EC_{50} = 0.8 \mu M$). TRAP-6 (0.3 and 0.6 mg/kg) has a triphasic effect on mean arterial blood pressure (MAP) in anesthetized rats with a short decrease, an increase, and then a longer decrease in MAP following intravenous administration.²

References

1. Vassallo, R.R., Jr., Kieber-Emmons, T., Chichowski, K., *et al.* Structure-function relationships in the activation of platelet thrombin receptors by receptor-derived peptides. *J. Biol. Chem.* **267**(9), 6081-6085 (1992).
2. Sicker, T., Wuchold, F., Kaiser, B., *et al.* Systemic vascular effects of thrombin and thrombin receptor activating peptide in rats. *Thromb. Res.* **101**(6), 467-475 (2001).

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, 12/20/2022

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