

PRODUCT INFORMATION



[Tyr⁰] Fibrinopeptide A (trifluoroacetate salt)

Item No. 44043

Formal Name:	N-L-tyrosyl-fibrinopeptide A (human), trifluoroacetate salt	
Synonym:	[Tyr ⁰] FPA	
Peptide Sequence:	YADSGEGDFLAEGGGVR-OH	H-Tyr-Ala-Asp-Ser-Gly-Glu-Gly-Asp-Phe-Leu-
MF:	C ₇₂ H ₁₀₆ N ₂₀ O ₂₈ • XCF ₃ COOH	Ala-Glu-Gly-Gly-Gly-Val-Arg-OH
FW:	1,699.8	
Purity:	≥95%	• XCF ₃ COOH
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

[Tyr⁰] Fibrinopeptide A (FPA) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the [Tyr⁰] FPA (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. [Tyr⁰] FPA (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

[Tyr⁰] FPA (trifluoroacetate salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

[Tyr⁰] Fibrinopeptide A (FPA) is a synthetic peptide derivative of FPA, which is generated via thrombin-induced cleavage of fibrinogen.¹ [Tyr⁰] FPA contains a tyrosine residue preceding the FPA sequence.

Reference

1. Coppens, M., Eikelboom, J.W., Gustafsson, D., *et al.* Translational success stories: Development of direct thrombin inhibitors. *Circ. Res.* **111(7)**, 920-929 (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, 09/26/2025

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