

PRODUCT INFORMATION



Fibrinogen (β chain) (41-62)-biotin Peptide

Item No. 39127

Overview and Properties

Contents:	This vial contains 1 mg of lyophilized peptide.
Synonyms:	Biotin-Phe-Ser-Ala-Arg-Gly-His-Arg-Pro-Leu-Asp-Lys-Lys-Arg-Glu-Glu-Ala-Pro-Ser-Leu-Arg-Pro-Ala, FGB
Peptide Sequence:	Biotin-FSARGHRPLDKKREEAPSLRPA
Uniprot No.:	P02675
Molecular Weight:	2,745.1
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥ 3 years

Description

Fibrinogen (β chain) (41-62)-biotin peptide is a biotinylated version of the secreted matrix protein fibrinogen, which undergoes citrullination and has been found as the citrullinated form in plasma circulating immune complexes of patients with rheumatoid arthritis.¹ Fibrinogen is subject to citrullination by peptidyl arginine deiminase 2 (PAD2) and PAD4, which can occur in the synovial fluid from patients with rheumatoid arthritis.^{2,3} Cayman's Fibrinogen (β Chain) (41-62)-biotin Peptide is intended for use as a negative control in the detection of autoantibodies against citrullinated fibrinogen and is the non-citrullinated form of Citrullinated Fibrinogen (R44 + R47 + R60) (41-62)-biotin Peptide (Item No. 39128)

References

1. Zhao, X., Okeke, N.L., Sharpe, O., *et al.* Circulating immune complexes contain citrullinated fibrinogen in rheumatoid arthritis. *Arthritis Res. Ther.* **10(4)**, R94 (2008).
2. Damiana, T., Damgaard, D., Sidemann, J.J., *et al.* Citrullination of fibrinogen by peptidylarginine deiminase 2 impairs fibrin clot structure. *Clin. Chim. Acta* **501**, 6-11 (2020).
3. Damgaard, D., Senolt, L., Nielsen, M.F., *et al.* Demonstration of extracellular peptidylarginine deiminase (PAD) activity in synovial fluid of patients with rheumatoid arthritis using a novel assay for citrullination of fibrinogen. *Arthritis Res. Ther.* **16(6)**, 498 (2014).

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, 12/14/2023

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