

PRODUCT INFORMATION

Citrullinated Fibrinogen (β chain) (R44 + R47 + R60) (41-62)-biotin Peptide Item No. 39128

Overview and Properties

Contents:	This vial contains 1 mg of lyophilized peptide.
Peptide Sequence:	Biotin-FSAXGHXPLDKKREEAPSLXPA (X=Citrulline)
Uniprot No.:	P02675
Molecular Weight:	2,747.9
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥ 3 years

Description

Citrullinated fibrinogen (β chain) (R44 + R47 + R60) (41-62)-biotin peptide is a biotinylated version of a tri-citrullinated form of the secreted matrix protein fibrinogen, which 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 Citrullinated Fibrinogen (β Chain) (R44 + R47 + R60) (41-62)-biotin Peptide is intended for the detection of autoantibodies against fibrinogen citrullinated at R44, R47, and R60.

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, 12/15/2023

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