

PRODUCT INFORMATION



CCP-biotin Peptide (non-cyclized)

Item No. 35876

Overview and Properties

Contents:	This vial contains contains 1 mg of lyophilized peptide.
Synonyms:	Citrullinated Filaggrin (Q447H, S449C, G459R, E460G, S462C) (R455) (447-467)-biotin Peptide (non-cyclized), Cyclic Citrullinated Peptide-biotin Peptide (non-cyclized)
Peptide Sequence:	Biotin-HQCHQESTXGRSRGRCSGS (X=Citrulline)
Molecular Weight:	2,569.7 kDa
Uniprot No.:	P20930
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥1 year

Description

CCP-biotin peptide (non-cyclized) is a biotinylated and non-cyclized version of circular citrullinated peptide (CCP).¹ The sequence corresponds to amino acids 447-467 of human filaggrin but contains Q447H, S449C, G459R, E460G, and S462C substitutions and is citrullinated at arginine 455. Cayman's CCP-biotin Peptide (non-cyclized) is intended for use in the detection of anti-citrulline autoantibodies, which may bind the linear peptide.

Reference

1. Won, P., Kim, Y., Jung, H., *et al.* Pathogenic role of circulating citrullinated antigens and anti-cyclic monoclonal citrullinated peptide antibodies in rheumatoid arthritis. *Front. Immunol.* **12**, 692242 (2021).

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, 11/28/2023

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

PRODUCT INFORMATION



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