

# PRODUCT INFORMATION



## CCP-biotin Peptide (non-cyclized)

Item No. 35876

### Overview and Properties

<b>Contents:</b>	This vial contains contains 1 mg of lyophilized peptide.
<b>Synonyms:</b>	Citrullinated Filaggrin (Q447H, S449C, G459R, E460G, S462C) (R455) (447-467)-biotin Peptide (non-cyclized), Cyclic Citrullinated Peptide-biotin Peptide (non-cyclized)
<b>Peptide Sequence:</b>	Biotin-HQCHQESTXGRSRGRCSRGS (X=Citrulline)
<b>Molecular Weight:</b>	2,569.7 Da
<b>Uniprot No.:</b>	P20930
<b>Form:</b>	Solid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Resuspension:</b>	Resuspend to 1 mg/ml in aqueous buffer of choice.

### Description

CCP-biotin peptide (non-cyclized) is a biotinylated and non-cyclized version of circular citrullinated peptide (CCP).<sup>1</sup> 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, 05/18/2026

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