

# PRODUCT INFORMATION



## NCP-biotin Peptide (non-cyclized)

Item No. 40255

### Overview and Properties

<b>Contents:</b>	This vial contains 1 mg of lyophilized peptide.
<b>Synonyms:</b>	Filaggrin (Q447H, S449C, G459R, E460G, S462C) (447-467)-biotin Peptide (non-cyclized), Non-citrullinated Peptide-biotin Peptide (non-cyclized)
<b>Peptide Sequence:</b>	Biotin-HQCHQESTRGRSRGRCGRSGS
<b>Molecular Weight:</b>	2,568.8 Da
<b>Uniprot No.:</b>	P20930
<b>Form:</b>	Lyophilized
<b>Storage:</b>	-20°C (as supplied); aliquot after resuspension and store at -20°C minimizing freeze/thaw
<b>Stability:</b>	≥1 year
<b>Resuspension:</b>	Resuspend to 1 mg/ml in water or aqueous buffer of choice.

### Description

NCP-biotin peptide (non-cyclized) is a biotinylated and non-cyclized version of non-citrullinated peptide (NCP).<sup>1</sup> The sequence corresponds to amino acids 447-467 of human filaggrin but contains Q447H, S449C, G459R, E460G, and S462C substitutions. Cayman's NCP-biotin Peptide (non-cyclized) is intended for use as a negative control and is the non-citrullinated version of CCP-biotin Peptide (non-cyclized) (Item No. 35876).

### 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, 05/07/2026

#### 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