

# PRODUCT INFORMATION



## CCP (trifluoroacetate salt)

Item No. 44486

**Formal Name:** ((4R,7S,13S,16S,19S,25S,28S,31S,34S,37S,40S,43R)-40-((1H-imidazol-4-yl)methyl)-43-((S)-5-amino-2-((S)-2-amino-3-(1H-imidazol-4-yl)propanamido)-5-oxopentanamido)-37-(3-amino-3-oxopropyl)-34-(2-carboxyethyl)-7,13,19-tris(3-guanidinopropyl)-28-((R)-1-hydroxyethyl)-16,31-bis(hydroxymethyl)-6,9,12,15,18,21,24,27,30,33,36,39,42-tridecaoxo-25-(3-ureidopropyl)-1,2-dithia-5,8,11,14,17,20,23,26,29,32,35,38,41-tridecaazacyclotetracontane-4-carbonyl)glycyl-L-arginyl-L-serylglycyl-L-seryl-L-lysine-2,2,2-trifluoroacetic acid

**Synonyms:** Citrullinated Filaggrin (Q447H, S449C, G459R, E460G, S462C, S468K) (R455) (447-468) (cyclized), Cyclic Citrullinated Peptide

**Peptide Sequence:** HQ[CHQESTXGRSRGRC]GRSGSK-OH  
(X=citrulline)

**MF:**  $C_{93}H_{157}N_{43}O_{33}S_2 \cdot XCF_3COOH$

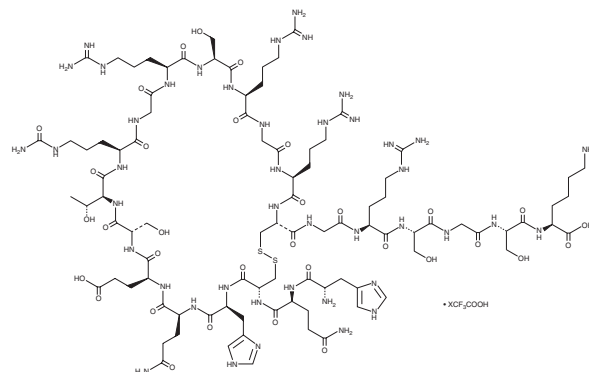
**FW:** 2,469.7

**Purity:**  $\geq 95$

**Supplied as:** A solid

**Storage:**  $-20^\circ C$

**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

CCP (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the CCP (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. CCP (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CCP (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. CCP (trifluoroacetate salt) is soluble ( $\geq 10$  mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

## Description

Cyclic citrullinated peptide (CCP) is a cyclic peptide that corresponds to amino acids 447-468 of human filaggrin but contains Q447H, S449C, G459R, E460G, S462C, and S468K substitutions and is citrullinated at arginine 455. This product can be used in the detection of anti-citrulline autoantibodies.

### 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, 12/02/2025

## 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