

PRODUCT INFORMATION

Tripeptide-10 Citrulline (trifluoroacetate salt)

Item No. 42215

Formal Name: L-lysyl-L- α -aspartyl-L-isoleucyl-N5-(aminocarbonyl)-L-ornithinamide, monotrifluoroacetate salt

Synonym: Lys-Asp-Ile-Cit-NH₂

Peptide Sequence: KDIX-NH₂ (X = citrulline)

MF: C₂₂H₄₂N₈O₇ • CF₃COOH

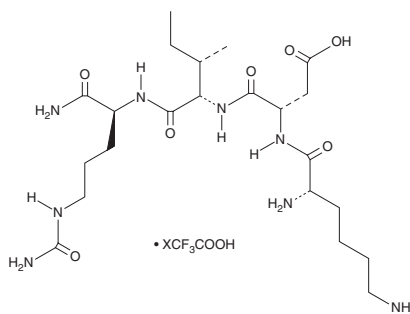
FW: 644.6

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Tripeptide-10 citrulline (trifluoroacetate salt) is supplied as a solid. Aqueous solutions of tripeptide-10 citrulline (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Tripeptide-10 citrulline (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Tripeptide-10 citrulline is a peptide inhibitor of collagen deformation.¹ It decreases collagen fibril diameter and fibrillogenesis in a synthetic human skin model when topically applied at a concentration of 0.01% (w/v). Formulations containing tripeptide-10 citrulline have been used in cosmetics.

Reference

1. Puig, A., Garcia Antón, J.M., and Mangues, M. A new decorin-like tetrapeptide for optimal organization of collagen fibres. *Int. J. Cosmet. Sci.* **30**(2), 97-104 (2008).

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, 01/17/2025

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