

PRODUCT INFORMATION

Collagen Type II (358-368)-biotin Peptide

Item No. 39125

Overview and Properties

Contents:	This vial contains 1 mg of lyophilized peptide.
Synonym:	Biotin-Gly-Ala-Arg-Gly-Leu-Thr-Gly-Arg-Pro-Gly-Asp-Ala
Peptide Sequence:	Biotin-GARGLTGRPGDA
Uniprot No.:	P02458
Molecular Weight:	1,353.5
Purity:	≥95%
Form:	A solid
Storage:	-20°C (as supplied)
Stability:	≥4 years
Resuspension:	Resuspend the peptide in water or buffer of choice at ≤1 mg/ml

Description

Collagen type II (358-368)-biotin peptide is a biotinylated version of the fibril-forming heterotrimeric protein collagen type II, which undergoes citrullination by peptidyl arginine deiminase 2 (PAD2) and PAD4.¹ Citrullinated collagen type II (358-368) has been found in synovial fluid isolated from patients with rheumatoid arthritis.^{1,2} Cayman's Collagen Type II (358-368)-biotin Peptide is intended for use as a negative control in the detection of autoantibodies against citrullinated collagen and is the non-citrullinated form of Citrullinated Collagen Type II (R360+R365) (358-368)-biotin Peptide (Item No. 39126).

References

1. Foulquier, C., Sebbag, M., Clavel, C., *et al.* Peptidyl arginine deiminase type 2 (PAD-2) and PAD-4 but not PAD-1, PAD-3, and PAD-6 are expressed in rheumatoid arthritis synovium in close association with tissue inflammation. *Arthritis Rheum.* **56(11)**, 3541-3553 (2007).
2. Uysal, H., Nandakumar, K.S., Kessel, C., *et al.* Antibodies to citrullinated proteins: Molecular interactions and arthritogenicity. *Immunol. Rev.* **233**, 9-33 (2009).

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, 10/14/2025

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