

PRODUCT INFORMATION

Citrullinated Collagen Type II (R360 + R365) (358-368)-biotin Peptide Item No. 39126

Overview and Properties

Synonym: Biotin-Gly-Ala-Arg(Cit)-Gly-Leu-Thr-Gly-Arg(Cit)-Pro-Gly-Asp-Ala
Peptide Sequence: Biotin-GAXGLTGXPGDA (X=Citrulline)
Purity: ≥95%
Supplied as: A solid
Storage: -20°C (as supplied)
Stability: ≥4 years

Reconstitution Instructions

Reconstitute in water or aqueous buffer of choice to 1 mg/ml. Upon reconstitution, store at 4°C for ≤1 month. For long term storage, aliquot and store at -20°C avoiding freeze/thaws.

Description

Citrullinated collagen type II (R360+R365) (358-368)-biotin peptide is a biotinylated derivative of a di-citrullinated peptide corresponding to a sequence found in the fibril-forming heterotrimeric protein collagen type II, which undergoes citrullination by peptidyl arginine deiminase 2 (PAD2) and PAD4.^{1,2} A synthetic peptide epitope derived from collagen type II citrullinated at R360 or R365 reacts with anti-citrullinated protein antibodies (ACPAs) isolated from mice with collagen-induced arthritis.² Citrullinated Collagen Type II (R360 + R365) (358-368)-biotin Peptide is intended for the detection of autoantibodies against collagen type II citrullinated at R360 and R365.

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**(1), 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, 11/20/2025

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