

PRODUCT INFORMATION



Citrullinated α -Enolase (R8 + R14) (1-19)-biotin Peptide

Item No. 37627

Overview and Properties

Contents:	This vial contains 1 mg of lyophilized peptide
Synonyms:	Citrullinated Enolase-1-biotin, α -Enolase Peptide (Citrulline Residues 8 + 14)
Peptide Sequence:	Biotin-SILKIHAXEIFDSXGNPTV (X=Citrulline)
Uniprot No.:	P06733
Molecular Weight:	2,381.5
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥3 years

Description

Citrullinated α -Enolase (R8 + R14) (1-19)-biotin peptide is a biotinylated and citrullinated form of a 19-amino acid N-terminal peptide fragment of α -enolase, a glycolytic enzyme that catalyzes the conversion of 2-phosphoglycerate to phosphoenolpyruvate.¹ α -Enolase also functions as a cell surface receptor for plasminogen on pathogens and activated immune cells, as an oxidative stress protein in endothelial cells, and as a chromatin binding partner to facilitate transcription.²⁻⁴ α -Enolase is an autoantigen in asthma, Hashimoto's encephalopathy, and rheumatoid arthritis, and has been found in the serum of pediatric patients with juvenile idiopathic arthritis.⁵⁻⁸ α -Enolase is also subject to citrullination by peptidyl arginine deiminases (PADs) and citrullinated α -enolase has been found in the synovial fluid of rheumatoid arthritis patients.⁷ Cayman's Citrullinated α -Enolase (R8 + R14) (1-19)-biotin Peptide is intended for use in the detection of autoantibodies against α -enolase citrullinated at R8 and R13.

References

1. Zhu, X., Miao, X., Wu, Y., *et al.* ENO1 promotes tumor proliferation and cell adhesion mediated drug resistance (CAM-DR) in non-Hodgkin's lymphomas. *Exp. Cell Res.* **335(2)**, 216-223 (2015).
2. Song, Y., Luo, Q., Long, H., *et al.* Alpha-enolase as a potential cancer prognostic marker promotes cell growth, migration, and invasion in glioma. *Mol. Cancer* **13**, 65 (2014).
3. Subramanian, A. and Miller, D.M. Structural analysis of α -enolase. Mapping the functional domains involved in down-regulation of the *c-myc* protooncogene. *J. Biol. Chem.* **275(8)**, 5958-5965 (2000).
4. Hsiao, K.-C., Shih, N.-Y., Fang, H.-L., *et al.* Surface α -enolase promotes extracellular matrix degradation and tumor metastasis and represents a new therapeutic target. *PLoS One* **8(7)**, e69354 (2013).
5. Nahm, D.-H., Lee, K.-H., Shin, J.-Y., *et al.* Identification of α -enolase as an autoantigen associated with severe asthma. *J. Allergy Clin. Immunol.* **118(2)**, 376-381 (2006).
6. Moore, T.L., Gillian, B.E., Crespo-Pagnussat, S., *et al.* Measurement and evaluation of isotypes of anti-citrullinated fibrinogen and anti-citrullinated α -enolase antibodies in juvenile idiopathic arthritis. *Clin. Exp. Rheumatol.* **32(5)**, 740-746 (2014).
7. Yoneda, M., Fujii, A., Ito, A., *et al.* High prevalence of serum autoantibodies against the amino terminal of α -enolase in Hashimoto's encephalopathy. *J. Neuroimmunol.* **185(1-2)**, 195-200 (2007).
8. Cong, Y., Wang, L., Peng, R., *et al.* Timosaponin AIII induces antiplatelet and antithrombotic activity via Gq-mediated signaling by the thromboxane A2 receptor. *Sci. Rep.* **6**, 38757 (2016).

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, 12/15/2023

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