

Rabbit Anti-Carbamyl-Lysine (CBL) Polyclonal Antibody

CATALOG NUMBER:	STA-078	STORAGE:	-20°C
QUANTITY AND CONCENTRATION:	50 µg of affinity purified antibody at 0.5 mg/mL in 75 mM PBS, pH 7.2, 75 mM NaCl, 0.5 mM EDTA and 0.02% NaN ₃		
SHELF LIFE:	1 year from date of receipt under proper storage conditions; aliquot to avoid multiple freeze thaw cycles		
HOST SPECIES:	Rabbit		
IMMUNOGEN:	CBL-KLH		
SPECIFICITY:	CBL-modified proteins		
APPLICATION:	Immunoblot (1:200 to 1:20,000) ELISA (1:200 to 1:20,000)		

Background

Carbamylation is a post-translational modification which occurs throughout the lifespan of proteins *in vivo*. Carbamylation results from the binding of isocyanic acid, spontaneously derived from high concentrations of urea and leading to the formation of carbamyl-lysine (CBL). The carbamylation of proteins is usually associated with a partial or complete loss of protein function. It is known that elevated urea directly induces the formation of potentially atherogenic carbamylated LDL (cLDL). High blood concentrations of urea leading to the carbamylation process were detected in uremic patients and patients with end-stage renal disease.

Recent Product Citations

1. Lee, Y.H. et al. (2020). Localisation of citrullinated and carbamylated proteins in inflamed gingival tissues from rheumatoid arthritis patients. *Clin Oral Investig.* doi: 10.1007/s00784-020-03452-9.
2. Lubbers, R. et al. (2019). Carbamylation reduces the capacity of IgG for hexamerisation and complement activation. *Clin Exp Immunol.* doi: 10.1111/cei.13411.
3. Verheul, M.K. et al. (2018). Pitfalls in the detection of citrullination and carbamylation. *Autoimmun Rev.* **17**(2):136-141. doi: 10.1016/j.autrev.2017.11.017.
4. Kluge-Beckerman, B. et al. (2016). Carbamylation of the amino-terminal residue (Gly1) of mouse serum amyloid A (SAA) promotes amyloid formation in a cell culture model. *FEBS Lett.* doi:10.1002/1873-3468.12472.
5. Skopelja, S. et al. (2016). The role for neutrophil extracellular traps in cystic fibrosis autoimmunity. *JCI Insight* **1**:e88912.
6. Challener, G. J. et al. (2016). Anti-carbamylated protein antibody levels correlate with anti-Sa (Citrullinated Vimentin) antibody levels in rheumatoid arthritis. *J Rheumatol.* **43**:273-81.

7. Bright, R. et al. (2015). Is there a link between carbamylation and citrullination in Periodontal Disease and Rheumatoid Arthritis?. *Med Hypotheses*. doi: 10.1016/j.mehy.2015.03.006.
8. Koro, C. et al. (2014). Carbamylation of immunoglobulin abrogates activation of the classical complement pathway. *Eur J Immunol*. **44**:3403-3412.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS 's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2013-2020: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.