

## CBL-BSA

**CATALOG NUMBER:** STA-379

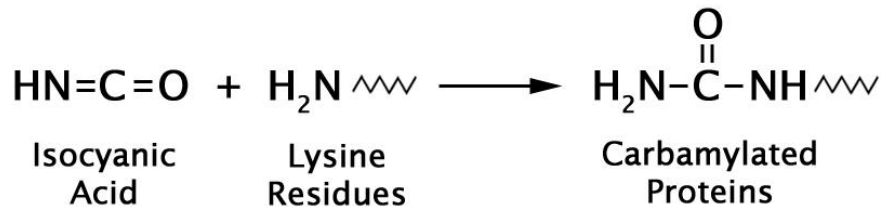
**STORAGE:** -20°C

**QUANTITY AND CONCENTRATION:** 100 µL of 0.1 mg/mL CBL-BSA in 1X PBS containing 0.05 mM EDTA, 0.001% NaN<sub>3</sub>, 5% Glycerol

**SHELF LIFE:** 1 year from date of receipt under proper storage conditions; aliquot to avoid multiple freeze thaw cycles

### 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) (Figure 1). 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.



**Figure 1: Formation of Carbamyl-Lysine (CBL) During Carbamylation of Proteins**

### Methods

Dilute the CBL-BSA with SDS-PAGE reducing sample buffer to 1.0-10 µg/mL and boil for 5 minutes. Load 10 µL per lane for western blot analysis of CBL protein adducts.

### References

1. Wang, Z., et al. (2007). *Nat. Med.* **13**:1176-1184.
2. Kraus, L.M., et al. (2001). *Swiss Med. Wkly.* **131**:139-144
3. Sirpal, S. (2009). *Clin. Sci. (Lond.)* **116**:681-695.
4. Balion, C.M., et al. (1998). *Kidney Int.* **53**:488-495.
5. Ascì, G., et al. (2008). *Nephrology (Carlton)* **13**:480-486.
6. Jaisson, S., et al. (2007). *FEBS Lett.* **581**:1509-1513.

7. Garnotel, R., et al. (2004). *FEBS Lett.* **563**:13-16.

### **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](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2012: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.