

# PRODUCT INFORMATION



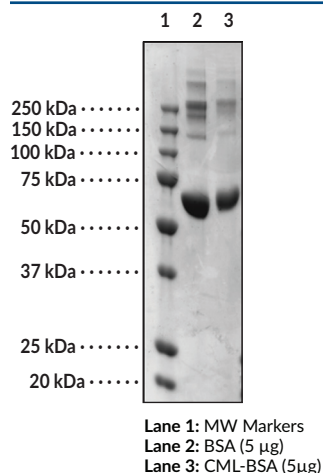
## CML-BSA

Item No. 22972

### Overview and Properties

<b>Synonyms:</b>	N <sup>ε</sup> -(1-Carboxymethyl)-L-lysine-Bovine Serum Albumin, N <sup>ε</sup> -(1-Carboxymethyl)-L-lysine-BSA
<b>Source:</b>	Albumin isolated from bovine plasma and modified by carboxymethylation
<b>Molecular Weight:</b>	69.3 kDa
<b>Storage:</b>	-20°C (as supplied) avoid freeze/thaw cycles by aliquoting the protein after resuspension
<b>Stability:</b>	≥2 years
<b>Supplied in:</b>	Lyophilized from a solution in PBS, pH 7.4, 200 µg per vial. Reconstitution with 0.2 ml of water will yield 1 mg/ml solution in 1X PBS.

### Image



### Description

Advanced glycation end products (AGEs) are formed from the nonenzymatic reaction of amino groups with reducing sugars.<sup>1-3</sup> AGEs have been implicated in diseases, such as diabetes mellitus, non-diabetic nephropathy, macrovascular disease, Alzheimer's disease, cataract, and ageing.<sup>1-3</sup> AGE receptors, such as the receptor for AGE (RAGE), mediate biological responses to AGEs, including endocytic uptake and degradation and induction of cytokines and growth factors.<sup>1-3</sup> CML-BSA was produced by incubating BSA with glyoxylic acid and sodium cyanoborohydride.

### References

1. Singh, R., Barden, A., Mori, T., et al. *Diabetologia* **44**(2), 129-46 (2001).
2. Giacco, F., and Brownlee, M. *Circ. Res.* **107**(9), 1058-1070 (2010).
3. Vistoli, G., De Maddis, D., Cipak, A., et al. *Free Radic. Res.* **47**(Suppl 1), 3-27 (2013).

**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, 09/20/2023

**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