

# PRODUCT INFORMATION



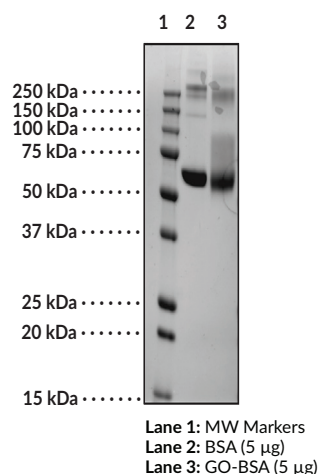
## GO-BSA

Item No. 22970

### Overview and Properties

**Synonyms:** Glyoxal-Bovine Serum Albumin, Glyoxal-BSA, GO-Bovine Serum Albumin  
**Source:** Albumin isolated from bovine plasma and modified with glyoxal  
**Molecular Weight:** 69.3 kDa  
**Storage:** -20°C (as supplied); avoid freeze/thaw cycles by aliquoting the protein after resuspension  
**Stability:** ≥2 years  
**Supplied in:** Lyophilized from a solution in PBS, pH 7.4, at 0.2 mg per vial. Reconstitution in 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> GO-BSA was produced by incubating BSA with glyoxal, followed by extensive dialysis.

### References

1. Singh, R., Barden, A., Mori, T., *et al.* Advanced glycation end-products: A review. *Diabetologia* **44**(2), 129-46 (2001).
2. Giacco, F. and Brownlee, M. Oxidative stress and diabetic complications. *Circ. Res.* **107**(9), 1058-1070 (2010).
3. Vistoli, G., De Maddis, D., Cipak, A., *et al.* Advanced glycoxidation and lipoxidation end products (AGEs and ALEs): an overview of their mechanisms of formation. *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