

PRODUCT INFORMATION



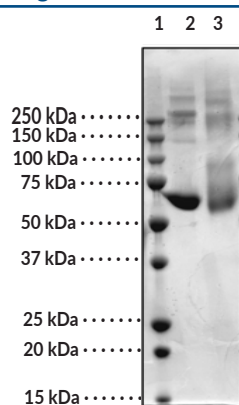
MG-BSA

Item No. 22971

Overview and Properties

Synonyms: Methylglyoxal-Bovine Serum Albumin, Methylglyoxal-BSA
Source: Albumin isolated from bovine plasma and modified with methylglyoxal
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
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Standards
Lane 2: BSA (5 µg)
Lane 3: MG-BSA (5 µg)

Representative gel image shown; actual purity may vary between each batch.

Description

Advanced glycation end products (AGEs) are formed from the nonenzymatic reaction of amino groups with reducing sugars.¹⁻³ AGEs have been implicated in diseases, such as diabetes mellitus, non-diabetic nephropathy, macrovascular disease, Alzheimer's disease, cataract, and ageing.¹⁻³ 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.¹⁻³ MG-BSA was produced by incubating BSA with methylglyoxal, 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, 08/25/2017

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