

PRODUCT INFORMATION



BSA-PEG(2000)

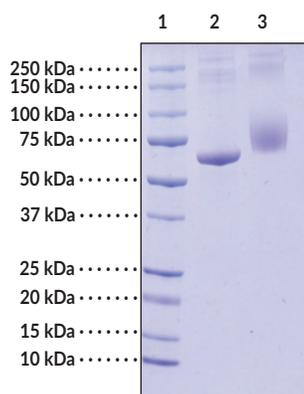
Item No. 43456

Overview and Properties

Synonyms: BSA-PEG(2K), m-Polyethylene Glycol-NHS ester BSA, PEGylated Bovine Serum Albumin
Source: Albumin isolated from bovine plasma and modified with methoxy-PEG(2000)-NHS ester
Molecular Weight: 66.5 kDa (unconjugated)
Storage: -20°C (as supplied)
Stability: ≥1 year
Supplied in: PBS, pH 7.4
Protein
Concentration: *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: BSA (4 μg)
Lane 3: BSA-PEG(2000) (4 μg)

SDS-PAGE Analysis of BSA-PEG(2000).

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

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, 03/18/2026

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

PRODUCT INFORMATION



Description

Polyethylene glycols (PEGs) are synthetic and hydrophilic polymers.^{1,2} They are linear or branched and contain a reactive end group, such as acrylate, methacrylate, dibenzocyclooctynol, or vinyl sulfonate, for covalent attachment to macromolecules or linkers. The opposite end group of PEGs is commonly a methyl group (methoxy PEG), however, hydroxy, amino, butoxy, and *tert*-butoxy end groups have also been used.¹ PEGs are non-toxic and are commonly used to prolong the *in vivo* circulation time of pharmaceutical agents, and PEGylated BSA has been used as a drug carrier *in vitro*.²⁻⁴ Free PEGs are non-immunogenic but become immunogenic when conjugated to a drug delivery nanosystem (DDS) or a macromolecule.¹ Immunogenicity of PEGs varies based on polymer length and branching, end group composition, and chemical nature of the PEG acceptor structure. Cayman's BSA-PEG(2000) protein is composed of methoxy-PEG(2000)-NHS ester and BSA at an approximately 10:1 molar ratio of methoxy-PEG(2000)-NHS ester:BSA. This product can be used as a positive control in ELISA and Western blot (WB) experiments using Cayman's Polyethylene Glycol Rabbit Monoclonal Antibody (Clone RM105) (Item No. 32180) and Polyethylene Glycol Rabbit Monoclonal Antibody - Biotinylated (Clone RM105) (Item No. 32381).

References

1. Kozma, G.T., Shimizu, T., Ishida, T., *et al.* Anti-PEG antibodies: Properties, formation, testing and role in adverse immune reactions to PEGylated nano-biopharmaceuticals. *Adv. Drug Deliv. Rev.* 154-155, 163-175 (2020).
2. Zhang, Z., Zhang, Y., Song, S., *et al.* Recent advances in the bioanalytical methods of polyethylene glycols and PEGylated pharmaceuticals. *J. Sep. Sci.* **43(9-10)**, 1978-1997 (2020).
3. Li, X.-Y., Li, T.-H., Guo, J.-S., *et al.* PEGylation of bovine serum albumin using click chemistry for the application as drug carriers. *Biotechnol. Prog.* **28(3)**, 856-861 (2012).
4. Ferrado, J.B., Perez, A.A., Menegon, M., *et al.* PEGylation of genistein-loaded bovine serum albumin nanoparticles and its effect on *in vitro* cell viability and genotoxicity properties. *Colloids Surf. B Biointerfaces* **222**, 113082 (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