

PRODUCT INFORMATION

BSA-Adrenate Polyunsaturated Fatty Acid Complex (1.67 mM)

Item No. 45953

Synonyms:	Bovine Serum Albumin-AdA, Bovine Serum Albumin-Adrenate, BSA-AdA, BSA-7(Z),10(Z),13(Z),16(Z)-Docosatetraenoic Acid
Supplied as:	1.67 mM Adrenate:0.8 mM BSA (2:1 adrenate:BSA) in 150 mM sodium chloride, pH 7.4
Storage:	-20°C
Stability:	≥2 years
Item Origin:	Animal/Bovine

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

Description

BSA-Adrenate Polyunsaturated Fatty Acid Complex (1.67 mM) is composed of adrenic acid (Item No. 90300) and fatty acid-free bovine serum albumin (BSA) at an approximately 2:1 molar ratio of adrenate:BSA. It was prepared under sterile conditions, then filtered and aliquoted into sterile vials. Cayman's BSA-Adrenate Polyunsaturated Fatty Acid Complex (1.67 mM) can be used for efficient fatty acid delivery to cells in culture for the purpose of monitoring lipid metabolism and inflammatory signaling pathways.^{1,2} It can also be used to sensitize cells to ferroptosis via its incorporation into phosphatidylethanolamines (PEs) and subsequent oxidation to form pro-ferroptotic PE species.³ It is suitable for use in short- and long-term cell culture applications (25+ hours). For best results, it is recommended that this product be used in conjunction with Cayman's BSA Control for BSA-Fatty Acid Complexes (5 mM) (Item No. 29556), prepared with fatty acid-free BSA.

References

1. Alsabeeh, N., Chausse, B., Kakimoto, P.A., *et al.* Cell culture models of fatty acid overload: Problems and solutions. *Biochim. Biophys. Acta Mol. Cell Biol. Lipids* **1863**(2), 143-151 (2018).
2. Brouwers, H., Jónasdóttir, H.S., Kuipers, M.E., *et al.* Anti-inflammatory and proresolving effects of the omega-6 polyunsaturated fatty acid adrenic acid. *J. Immunol.* **205**(10), 2840-2849 (2020).
3. Kagan, V.E., Mao, G., Qu, F., *et al.* Oxidized arachidonic and adrenic PEs navigate cells to ferroptosis. *Nat. Chem. Biol.* **13**(1), 81-90 (2017).

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, 06/30/2026

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