

# PRODUCT INFORMATION

## BSA-Eicosapentaenoate Polyunsaturated Fatty Acid Complex (1 mM)

Item No. 43473

### Overview and Properties

<b>Synonyms:</b>	Bovine Serum Albumin-Eicosapentaenoate, Bovine Serum Albumin-EPA, BSA-EPA
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Supplied in:</b>	1 mM EPA:0.17 mM BSA in 150 mM sodium chloride
<b>Item Origin:</b>	Animal/Bovine

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

### Description

BSA-Eicosapentaenoate polyunsaturated fatty acid complex is composed of eicosapentaenoic acid (EPA; Item Nos. 90110 | 90110.1) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of eicosapentaenoate:BSA. It has been used for efficient fatty acid delivery to cells in culture for the purpose of monitoring fatty acid oxidation or similar processes in various cellular metabolic studies.<sup>1-3</sup> BSA/BSA-Fatty acids are suitable for use in short-term cell culture applications (acute treatment to 18 hours); however, for long-term applications (25+ hours) the product should be filter-sterilized using a 0.2 µm filter and sterile receptacle, which will not affect its performance. For best results, it is recommended that this product be used in conjunction with BSA control for BSA-fatty acid complexes (1 mM) (Item No. 34932).

### 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. Wiesenfeld, P.W., Babu, U.S., and O'Donnell, M.W. Effect of long-chain fatty acids in the culture medium on fatty acid composition of WEHI-3 and J774A.1 cells. *Comp. Biochem. Physiol. B Biochem. Mol. Biol.* **128**(1), 123-134 (2001).
3. Arita, K., Kobuchi, H., Utsumi, T., *et al.* Mechanism of apoptosis in HL-60 cells induced by n-3 and n-6 polyunsaturated fatty acids. *Biochem. Pharmacol.* **62**(7), 821-828 (2001).

**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, 11/21/2025

**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