

# PRODUCT INFORMATION



## BSA-Linoleate Reagent Set (1 mM)

Item No. 44717

### Overview and Properties

**Synonyms:** Bovine Serum Albumin-LA, Bovine Serum Albumin-Linoleate, BSA-LA  
**Supplied as:** Item No. 44718: 1 mM linoleate:0.17 mM BSA in 150 mM sodium chloride, pH 7.4  
Item No. 44719: 0.17 mM BSA in 150 mM sodium chloride, pH 7.4  
**Sterility:** Sterile

### Materials Supplied

Item Number	Item Name	Quantity/Size	Storage Temperature
44718	BSA-Linoleate Complex (1 mM)	1 vial/5 ml	-20°C
44719	BSA Control (Linoleate Complex)	1 vial/5 ml	-20°C

### Storage and Stability

Store individual components as directed in the Materials Supplied section and use before the expiration date indicated on the label.

### Description

The BSA-Linoleate Reagent Set contains a 1 mM solution of BSA-linoleate complex and a fatty acid-free BSA control. BSA-Linoleate Reagent Set (1 mM) is composed of linoleic acid (Item Nos. 90150 | 90150.1 | 21909) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of linoleate:BSA. The BSA-linoleate complex and BSA control were prepared at the same time with the same lot of BSA under sterile conditions, then filtered and aliquoted into sterile vials. BSA-linoleate complexes have 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>1-4</sup> Cayman's BSA-Linoleate Reagent Set is suitable for use in short- and long-term cell culture applications (25+ hours).

### 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. Menon, D., Salloum, D., Bernfeld, E., *et al.* Lipid sensing by mTOR complexes via *de novo* synthesis of phosphatidic acid. *J. Biol. Chem.* **292**(15), 6303-6311 (2017).
3. 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).
4. Colquhoun, A. and Curi, R. Regulation of tumour cell fatty acid oxidation by *n-6* polyunsaturated fatty acids. *Biochem. Soc. Trans.* **25**(4), S681 (1997).

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, 12/17/2025

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