

PRODUCT INFORMATION



BSA-Oleate Reagent Set (5 mM)

Item No. 44712

Overview and Properties

Synonyms: Bovine Serum Albumin-OA, Bovine Serum Albumin-Oleate, BSA-OA
Supplied as: Item No. 44713: 5 mM Oleate:0.8mMBSA(6:1 oleate:BSA) in 150 mM sodium chloride, pH 7.4
Item No. 44714: 0.8 mM BSA in 150 mM sodium chloride, pH 7.4
Sterility: Sterile

Materials Supplied

Item Number	Item Name	Quantity/Size	Storage Temperature
44713	BSA-Oleate Complex (5 mM)	1 vial/5 ml	-20°C
44714	BSA Control (Oleate 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-Oleate Reagent Set contains a 5 mM solution of BSA-oleate complex and a fatty acid-free BSA control. BSA-Oleate Complex (5 mM) is composed of oleic acid (Item Nos. 90260 | 24659) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of oleate:BSA. The BSA-oleate 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-oleate 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.¹⁻³ Cayman's BSA-Oleate Reagent Set is suitable for use in short- and long-term cell culture applications (25+ hours).

References

- 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).
- Trigatti, B.L. and Gerber, G.E. A direct role for serum albumin in the cellular uptake of long-chain fatty acids. *Biochem. J.* **308**(Pt. 1), 155-159 (1995).
- Chausse, B., Kakimoto, P.A., Caldeira-da-Silva, C.C., *et al.* Distinct metabolic patterns during microglial remodeling by oleate and palmitate. *Biosci. Rep.* **39**(4), BSR20190072 (2019).

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, 12/16/2025

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