

PRODUCT INFORMATION



BSA-Stearate Saturated Fatty Acid Complex (1 mM)

Item No. 38107

Synonym: Bovine Serum Albumin-SA, Bovine Serum Albumin-Stearate, BSA-Octadecanoic Acid, BSA-SA

Supplied as: 1 mM Stearate:0.17 mM BSA in 150 mM sodium chloride

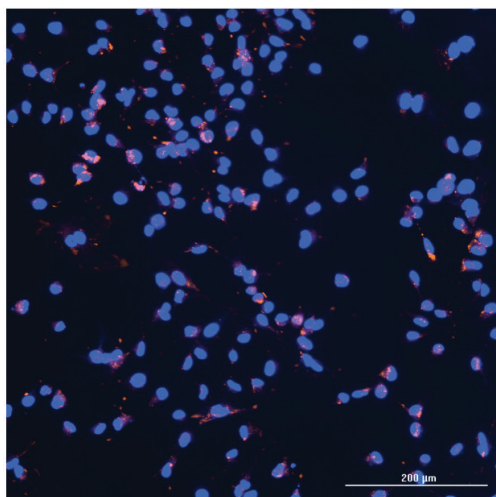
Storage: -20°C (as supplied)

Stability: ≥2 years

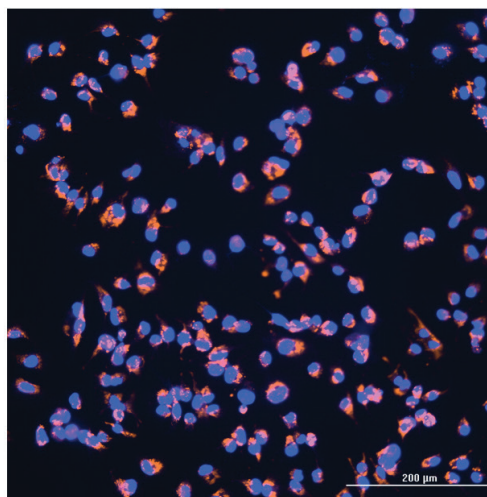
Item Origin: Animal/Bovine

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

Image



BSA Control



BSA-Stearate

Huh7 hepatocytes were plated and treated with 200 μ M BSA-Stearate Saturated Fatty Acid Complex (Item No. 38107) or BSA Control for BSA-Fatty Acid Complexes (Item No. 34932) for 24 hours. On day two, cells were incubated with Hoechst (blue) and Nile Red (red) for 15 minutes before imaging, showing lipid droplet accumulation.

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, 07/13/2026

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

BSA-Stearate Saturated Fatty Acid Complex (1 mM) is composed of stearic acid (Item No. 10011298) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of stearate:BSA. It was prepared under sterile conditions, then filtered and aliquoted into sterile vials. Cayman's BSA-Stearate Saturated Fatty Acid Complex (1 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 BSA-stearate complexes can also be used in colony formation, migration, and senescence assays to study the role of stearic acid in cancer *in vitro*.² Cayman's BSA-Stearate Saturated Fatty Acid Complex (1 mM) 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 (1 mM) (Item No. 34932), 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. Chen, P., Tian, J., Zhou, Y., *et al.* Metabolic flux analysis reveals the roles of stearate and oleate on CPT1C-mediated tumor cell senescence. *Int. J. Biol. Sci.* **19**(7), 2067-2080 (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