

PRODUCT INFORMATION



BSA-Octanoate Saturated Fatty Acid Complex (1 mM)

Item No. 34933

Overview and Properties

Storage: -20°C (as supplied)
Stability: ≥1 year
Supplied in: 1 mM octanoate : 0.17 mM BSA in 150 mM sodium chloride, pH 7.4
Item Origin: Animal/Bovine

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

Description

BSA-Octanoate saturated fatty acid complex is composed of octanoic acid (Item No. 33674) and bovine serum albumin (BSA) at an approximately 6:1 molar ratio of octanoate:BSA. BSA-Octanoate complexes can be 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.^{1,2} The BSA/BSA-FAs are acceptable for use with short-term cell culture applications (acute treatment to 18 hours), however, for long-term applications (25+ hours) the product can/should be filter sterilized using a 0.2 µm filter and sterile receptacle without affecting the 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. Tachibana, S., Sato, K., Cho, Y., *et al.* Octanoate reduces very low-density lipoprotein secretion by decreasing the synthesis of apolipoprotein B in primary cultures of chicken hepatocytes. *Biochim. Biophys. Acta* **1737**(1), 36-43 (2005).
2. Guo, W., Xie, W., and Han, J. Modulation of adipocyte lipogenesis by octanoate: Involvement of reactive oxygen species. *Nutr. Metab. (Lond.)* **3**, 30 (2006).

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, 05/10/2023

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