

PRODUCT INFORMATION



BSA Control for BSA-Fatty Acid Complexes (1 mM)

Item No. 34932

Supplied as: 0.17 mM BSA in 150 mM sodium chloride, pH 7.4
Storage: -20°C; avoid freeze/thaw cycles by aliquoting protein
Stability: ≥1 year
Item Origin: Animal/Bovine

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

Description

BSA control for BSA-fatty acid complexes (1 mM) is a fatty acid-free preparation of BSA intended for use as a control for BSA-fatty acid complexes containing 1 mM fatty acid (Item Nos. 34931 | 34933). It can be used as a control for experiments using BSA-fatty acid complexes to deliver fatty acids to cells in culture for the purpose of monitoring fatty acid oxidation or similar processes in various cellular metabolic studies.¹⁻³ 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.

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. Wang, D., Green, M.F., McDonnell, E., *et al.* Oxygen flux analysis to understand the biological function of sirtuins. *Methods Mol. Biol.* **1077**, 241-258 (2013).
3. Pardo, V., González-Rodríguez, Á., Guijas, C., *et al.* Opposite cross-talk by oleate and palmitate on insulin signaling in hepatocytes through macrophage activation. *J. Biol. Chem.* **290(18)**, 11663-11677 (2015).

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, 09/13/2021

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