

PRODUCT INFORMATION

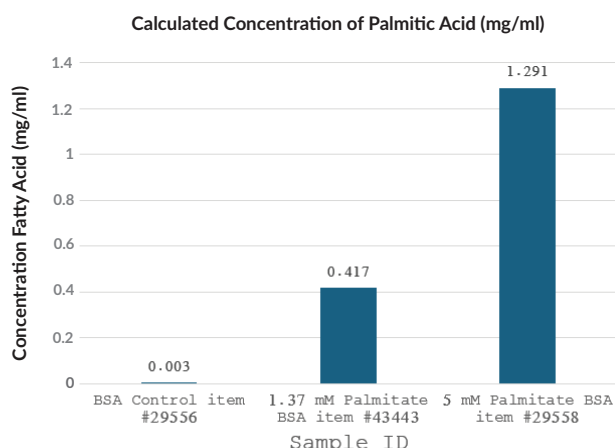
BSA-Palmitate Saturated Fatty Acid Complex (5 mM)

Item No. 29558

Synonym: Bovine Serum Albumin-PA, Bovine Serum Albumin-Palmitate, BSA-PA
Supplied as: 5 mM Palmitate:0.8 mM BSA (6:1 palmitate:BSA) in 150 mM sodium chloride, pH 7.4
Storage: -20°C
Stability: ≥1 year
Item Origin: Animal/Bovine

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

Images



Fatty acid concentration in BSA-Palmitate Saturated Fatty Acid Complex (1.67 mM) (Item No. 43443) and BSA-Palmitate Saturated Fatty Acid Complex (5 mM) (Item No. 29558) compared to BSA Control for BSA-Fatty Acid Complexes (5 mM) (Item No. 29556) using GC-MS analysis.

Description

BSA-Palmitate saturated fatty acid complex is composed of palmitic acid (Item No. 10006627) and bovine serum albumin (BSA) at an approximately 6:1 molar ratio of palmitate:BSA. It 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.¹⁻³ 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. For best results, it is recommended that this product be used in conjunction with BSA control for fatty acid complexes (5 mM) (Item No. 29556).

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).
- 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).
- Bentebibel, A., Sebastián, D., Herrero, L., et al. Novel effect of C75 on carnitine palmitoyltransferase I activity and palmitate oxidation. *Biochemistry* **45**(14), 4339-4350 (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, 06/03/2025

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