

# PRODUCT INFORMATION



## BSA-Palmitate Reagent Set (5 mM)

Item No. 44698

### Overview and Properties

**Synonyms:** Bovine Serum Albumin-PA, Bovine Serum Albumin-Palmitate, BSA-PA  
**Supplied as:** Item No. 44699: 5 mM Palmitate:0.8 mM BSA (6:1 palmitate:BSA) in 150 mM sodium chloride, pH 7.4  
Item No. 44670: 0.8 mM BSA in 150 mM sodium chloride, pH 7.4  
**Sterility:** Sterile

### Materials Supplied

| Item Number | Item Name                       | Quantity/Size | Storage Temperature |
|-------------|---------------------------------|---------------|---------------------|
| 44699       | BSA-Palmitate Complex (5 mM)    | 1 vial/5 ml   | -20°C               |
| 44700       | BSA Control (Palmitate 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-Palmitate Reagent Set contains a 5 mM solution of BSA-palmitate complex and a fatty acid-free BSA control. BSA-Palmitate Complex (5 mM) is composed of palmitic acid (Item No. 10006627) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of palmitate:BSA. The BSA-palmitate 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-palmitate 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.<sup>1-3</sup> Cayman's BSA-Palmitate 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).
- 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, 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