

# PRODUCT INFORMATION



## HSA-Palmitate Saturated Fatty Acid Complex (5 mM)

Item No. 45018

<b>Synonym:</b>	Human Serum Albumin-PA, Human Serum Albumin-Palmitate, HSA-PA
<b>Supplied as:</b>	5 mM Palmitate:0.8 mM HSA (6:1 palmitate:HSA) in 150 mM sodium chloride, pH 7.4
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥2 years
<b>Item Origin:</b>	Human

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

### Description

HSA-Palmitate saturated fatty acid complex is composed of palmitic acid (Item No. 10006627) and fatty acid-free human serum albumin (HSA) at an approximately 6:1 molar ratio of palmitate:HSA. It has 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> HSA/HSA-Fatty acids are suitable for use in short-term cell culture applications (acute treatment to 18 hours); however, for long-term applications (25+ hours) the product should be filter-sterilized using a 0.2 µm filter and sterile receptacle, which will not affect its performance. For best results, it is recommended that this product be used in conjunction with HSA control for HSA-fatty acid complexes (5 mM) (Item No. 45019).

### References

1. Spector, A.A., John, K., and Fletcher, J.E. Binding of long-chain fatty acids to bovine serum albumin. *J. Lipid Res.* **10(1)**, 56-67 (1969).
2. Twei, V.C., Ha, J.-S., and Ha, C.-E. Effects of human serum albumin complexed with free fatty acids on cell viability and insulin secretion in the hamster pancreatic β-cell line HIT-T15. *Life Sci.* **88(17-18)**, 810-818 (2011).
3. Simard, J.R., Zunszain, P.A., Ha, C.-E., et al. Locating high-affinity fatty acid-binding sites on albumin by x-ray crystallography and NMR spectroscopy. *Proc. Natl. Acad. Sci. U.S.A.* **102(50)**, 17958-17963 (2005).

#### 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, 01/22/2026

#### 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