

PRODUCT INFORMATION

HSA Control for HSA-Fatty Acid Complexes (5 mM)

Item No. 45019

Overview and Properties

Synonyms:	Human Serum Albumin
Supplied in:	0.8 mM HSA in 150 mM sodium chloride, pH 7.4
Storage:	-20°C (as supplied)
Stability:	≥2 years
Item Origin:	Human

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

Description

HSA control for HSA-fatty acid complexes (5 mM) is a fatty acid-free preparation of human serum albumin (HSA) intended for use as a control for HSA-fatty acid complexes containing 5 mM fatty acid (Item No. 45018). It can be used as a control for experiments using HSA-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 The HSA/HSA-FAs 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.

Reference

1. Tuei, 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).

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, 01/23/2026

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