

PRODUCT INFORMATION

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

Item No. 43649

Overview and Properties

Synonym: Human Serum Albumin
Source: Human
Molecular Weight: 66.5 kDa
Storage: -20°C (as supplied)
Stability: ≥2 years
Supplied in: 0.17 mM HSA in 150 mM sodium chloride, pH 7.4

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

Description

HSA control for HSA-fatty acid complexes (1 mM) is a fatty acid-free preparation of human serum albumin (HSA) intended for use as a control for HSA-fatty acid complexes containing 1 mM fatty acid (Item No. 43472). 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