

# PRODUCT INFORMATION

## BSA-13(Z)-Docosenoate Monounsaturated Fatty Acid Complex (1 mM) Item No. 39150

<b>Synonym:</b>	BSA-Erucate
<b>Supplied as:</b>	1 mM Erucate:0.17 mM BSA in 150 mM sodium chloride, pH 7.4
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥1 year
<b>Item Origin:</b>	Animal/Bovine

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

### Description

---

BSA-13(Z)-docosenoate monounsaturated fatty acid (FA) complex is composed of 13(Z)-docosenoic acid (Item Nos. 90175 | 38975) and bovine serum albumin (BSA) at an approximately 6:1 molar ratio of erucate:BSA. 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 BSA-fatty acid complexes (1 mM) (Item No. 34932).

#### 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, 10/18/2023

#### 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](http://WWW.CAYMANCHEM.COM)