

PRODUCT INFORMATION

BSA-cis-Parinarate Polyunsaturated Fatty Acid Complex (1 mM)

Item No. 44413

Synonyms:	Bovine Serum Albumin-cis-Parinarate, BSA-9(Z),11(E),13(E),15(Z)-Octadecatetraenoic Acid
Ex./Em. Max:	320/432 nm
Supplied as:	1 mM cis-parinarate:0.17 mM BSA (6:1 cis-parinarate:BSA) in 150 mM sodium chloride, pH 7.4
Storage:	-20°C
Stability:	≥2 years
Item Origin:	Animal/Bovine

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

Description

BSA-cis-Parinarate Polyunsaturated Fatty Acid Complex (1 mM) is composed of cis-parinaric acid (Item No. 71430) and fatty acid-free bovine serum albumin (BSA) at an approximately 6:1 molar ratio of cis-parinarate:BSA. It can be used for efficient fatty acid delivery to cells in culture for the purpose of monitoring lipid metabolism and inflammatory signaling pathways.¹⁻³ In cell culture media, cis-parinaric acid displays excitation/emission maxima of 320/432 nm, respectively, and can be used to measure phospholipase and lipase activity or to monitor lipid peroxidation.⁴⁻⁸ Cayman's BSA-cis-Parinarate Polyunsaturated Fatty Acid Complex (1 mM) is 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 Cayman's BSA Control for BSA-Fatty Acid Complexes (1 mM) (Item No. 34932), prepared with fatty acid-free BSA.

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. Tsei, 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).
4. Gomes, A., Fernandes, E., and Lima, J.L.F.C. Fluorescence probes used for detection of reactive oxygen species. *J. Biochem. Biophys. Methods* **65**(2-3), 45-80 (2005).
5. Wolf, C., Sagaert, L., and Berezat, G. A sensitive assay of phospholipase using the fluorescent probe 2-parinaroyllecithin. *Biochem. Biophys. Res. Commun.* **99**(1), 275-283 (1981).
6. Beisson, F., Ferté, N., Nari, J., et al. Use of naturally fluorescent triacylglycerols from *Parinari glaberrimum* to detect low lipase activities from *Arabidopsis thaliana* seedlings. *J. Lipid Res.* **40**(12), 2313-2321 (1999).
7. McGuire, S.O., James-Kracke, M.R., Sun, G.Y., et al. An esterification protocol for cis-parinaric acid-determined lipid peroxidation in immune cells. *Lipids* **32**(2), 219-226 (1997).
8. de Hingh, Y.C.M., Meyer, J., Fischer, J.C., et al. Direct measurement of lipid peroxidation in submitochondrial particles. *Biochemistry* **34**(39), 12755-12760 (1995).

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, 06/16/2026

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