

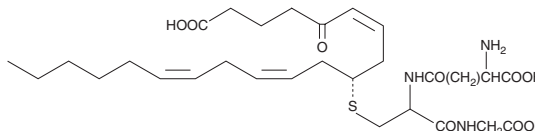
PRODUCT INFORMATION



FOG9

Item No. 34265

Formal Name: 5-oxo-6Z,9S-(S-glutathionyl)-11Z,14Z-eicosatrienoic acid
Synonym: 5-OxoETE-glutathione
MF: C₃₀H₄₇N₃O₉S
FW: 625.8
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 276 nm
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥2 years



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

Laboratory Procedures

FOG9 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of FOG9 in these solvents is approximately 50 mg/ml.

FOG9 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of FOG9 should be diluted with the aqueous buffer of choice. The solubility of FOG9 in PBS (pH 7.2) is approximately 100 µg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

FOG9 is formed through the conjugation of 5-oxoETE with glutathione in a 1,6 Michael addition at C-9. The biological activity of FOG9 has not been determined. However, FOG7, a 1,4-adduct formed within murine peritoneal macrophages, was found to be highly potent in stimulating eosinophil and neutrophil chemotaxis.¹ FOG7 was also capable of initiating actin polymerization without elevating intracellular free calcium ion concentration within either the eosinophil or polymorphonuclear.¹

Reference

1. Bowers, R.C., Hevko, J., Henson, P.M., et al. A novel glutathione containing eicosanoid (FOG7) chemotactic for human granulocytes. *J. Biol. Chem.* 275, 29931-29934 (2000).

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, 03/07/2024

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