

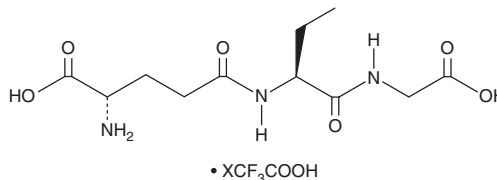
PRODUCT INFORMATION



Ophthalmic Acid (trifluoroacetate salt)

Item No. 35343

Formal Name: L-γ-glutamyl-(2S)-2-aminobutanoyl-glycine, trifluoroacetate salt
MF: C₁₁H₁₉N₃O₆ • XCF₃COOH
FW: 289.3
Purity: ≥98%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Ophthalmic acid (trifluoroacetate salt) 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 ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of ophthalmic acid (trifluoroacetate salt) in ethanol is approximately 12 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Ophthalmic acid (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of ophthalmic acid (trifluoroacetate salt) should be diluted with the aqueous buffer of choice. The solubility of ophthalmic acid (trifluoroacetate salt) in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ophthalmic acid is a glutathione (GSH) analog and biomarker of hepatic glutathione depletion.¹ It is formed *via* the GSH synthesis pathway, substituting 2-aminobutyrate for cysteine. Plasma and hepatic levels of ophthalmic acid are increased in fasted mice and in a mouse model of acetaminophen-induced hepatotoxicity.^{1,2}

References

1. Kobayashi, S., Lee, J., Takao, T., *et al.* Increased ophthalmic acid production is supported by amino acid catabolism under fasting conditions in mice. *Biochem. Biophys. Res. Commun.* **491(3)**, 649-655 (2017).
2. Soga, T., Baran, R., Suematsu, M., *et al.* Differential metabolomics reveals ophthalmic acid as an oxidative stress biomarker indicating hepatic glutathione consumption. *J. Biol. Chem.* **281(24)**, 16768-16776 (2006).

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, 12/01/2025

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