

PRODUCT INFORMATION



Papain Inhibitor (trifluoroacetate salt)

Item No. 31404

Formal Name: glycyglycyl-L-tyrosyl-L-arginine, trifluoroacetate salt

Synonyms: GGYR, Gly-Gly-Tyr-Arg

MF: C₁₉H₂₉N₇O₆ • CF₃COOH

FW: 565.5

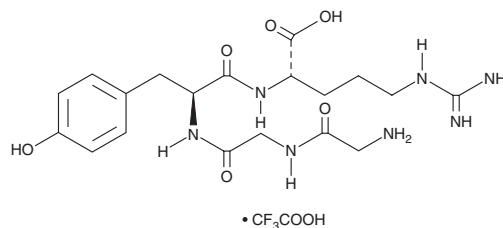
Purity: ≥95%

UV/Vis.: λ_{max}: 227 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Papain inhibitor (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the papain inhibitor (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Papain inhibitor (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of papain inhibitor (trifluoroacetate salt) in these solvents is approximately 10 mg/ml in ethanol and 1 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of papain inhibitor (trifluoroacetate salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of papain inhibitor in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Papain inhibitor is a synthetic peptide inhibitor of the cysteine protease papain ($K_i = 113 \mu\text{M}$).¹

Reference

1. Funk, M.O., Nakagawa, Y., Skochopole, J., et al. Affinity chromatographic purification of papain. *Int. J. Pept. Protein Res.* **13**(3), 296-303 (1979).

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, 12/12/2022

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