

PRODUCT INFORMATION



Mastoparan X (trifluoroacetate salt)

Item No. 36903

CAS Registry No.: 72093-22-2

Formal Name: L-isoleucyl-L-asparaginyl-L-tryptophyl-L-lysylglycyl-L-isoleucyl-L-alanyl-L-alanyl-L-methionyl-L-alanyl-L-lysyl-L-lysyl-L-leucyl-L-leucinamide, trifluoroacetate salt

Synonyms: Mastoparan V, MPV, MPX

Peptide Sequence: INWKGIAMAKKLL-NH₂

MF: C₇₃H₁₂₆N₂₀O₁₅S • XCF₃COOH

FW: 1,556.0

Purity: ≥98%

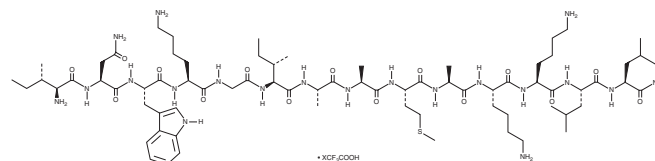
UV/Vis.: λ_{max}: 221 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Synthetic



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

Laboratory Procedures

Mastoparan X (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the mastoparan X (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Mastoparan X (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of mastoparan X (trifluoroacetate salt) in ethanol is approximately 50 mg/ml and approximately 25 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 mastoparan X (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of mastoparan X (trifluoroacetate salt) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Mastoparan X is an antimicrobial peptide that has been found in wasp venom.¹ It is active against several Gram-positive and Gram-negative bacteria (MICs = 3-64 and 4-256 µg/ml, respectively) and induces permeabilization of *E. coli* cytoplasmic membranes by 86% when used at a concentration of 6.4 µg/ml. Mastoparan X induces GDP/GTP exchange by G_i in a concentration-dependent manner and binds to calmodulin (K_d = ~0.9 nM) in cell-free assays.^{2,3} It induces necrosis in T98G glioblastoma multiforme (GBM) cells when used at a concentration of 20 µM.⁴ Mastoparan X (128 µg/ml) induces degranulation of isolated rat peritoneal mast cells and hemolysis of isolated sheep, chicken, and human erythrocytes in a concentration-dependent manner.¹ *In vivo*, mastoparan X (20 mg/kg) increases survival, decreases disease severity, and prevents damage to jejunal villi and crypts in a mouse model of enterohemorrhagic *E. coli* infection.⁵

References

1. Lin, C., -H., Tzen, J.T.C., Shyu, C.-L., et al. *Peptides* **32**(10), 2027-2036 (2011).
2. Malencik, D.A. and Anderson, S.R. *Biochem. Biophys. Res. Commun.* **114**(1), 50-56 (1983).
3. Higashijima, T., Burnier, J., and Ross, E.M. *J. Biol. Chem.* **265**(24), 14176-14186 (1990).
4. da Silva, A.M., Silva-Gonçalves, L.C., Oliveira, F.A., et al. *Mol. Neurobiol.* **55**(7), 5490-5504 (2018).
5. Zhao, X., Wang, L., Zhi, C., et al. *Front. Microbiol.* **12**, 644887 (2021).

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, 08/03/2023

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