

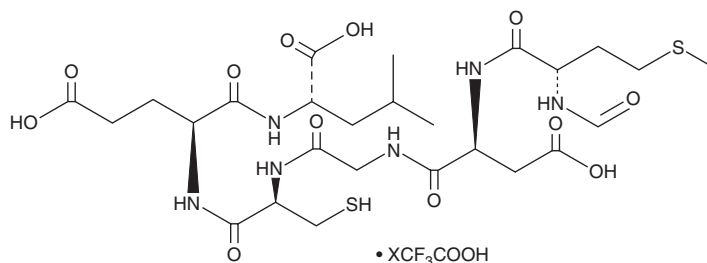
PRODUCT INFORMATION



Foxy-5 (trifluoroacetate salt)

Item No. 30983

Formal Name: N-formyl-L-methionyl-L- α -aspartylglycyl-L-cysteinyl-L- α -glutamyl-L-leucine
Synonyms: N-formyl-Met-Asp-Gly-Cys-Glu-Leu, N-formyl-MDGCEL
MF: C₂₆H₄₂N₆O₁₂S₂ • XCF₃COOH
FW: 694.8
Purity: ≥98%
Supplied as: A 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

Foxy-5 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the foxy-5 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Foxy-5 (trifluoroacetate salt) is soluble in ethanol.

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 foxy-5 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of foxy-5 (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

Foxy-5 is a Wnt5a peptide mimetic with anticancer activity.^{1,2} It increases adhesion and reduces cell motility of MDA-MB-468 breast cancer cells when used at concentrations ranging from 1 to 1,000 μ M.¹ Foxy-5 (2 mg/kg) reduces the number of lymph node metastases without affecting primary tumor growth in a DU145-Luc mouse xenograft model.²

References

1. Säfholm, A., Leandersson, K., Dejmek, J., *et al.* A formylated hexapeptide ligand mimics the ability of Wnt-5a to impair migration of human breast epithelial cells. *J. Biol. Chem.* **281**(5), 2740-2749 (2006).
2. Canesin, G., Evans-Axelsson, S., Hellsten, R., *et al.* Treatment with the WNT5A-mimicking peptide Foxy-5 effectively reduces the metastatic spread of WNT5A-low prostate cancer cells in an orthotopic mouse model. *PLoS One* **12**(9), e0184418 (2017).

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, 11/02/2022

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