

PRODUCT INFORMATION



cyclo(RGDyC) (trifluoroacetate salt)

Item No. 34945

Formal Name: 2-((2S,5R,8R,11S)-11-(3-guanidinopropyl)-5-(4-hydroxybenzyl)-8-(mercaptomethyl)-3,6,9,12,15-pentaoxo-1,4,7,10,13-pentaazacyclopentadecan-2-yl)acetic acid, trifluoroacetate salt

Synonyms: c(RGDyC), cyclo(Arg-Gly-Asp-D-Tyr-Cys)

MF: C₂₄H₃₄N₈O₈S • XCF₃COOH

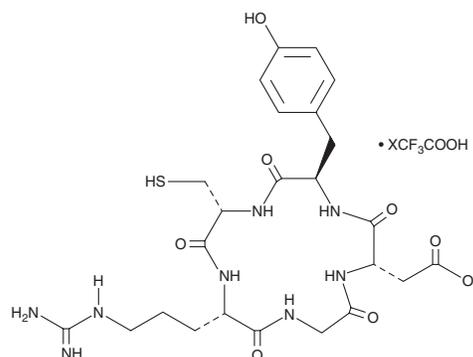
FW: 594.6

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

cyclo(RGDyC) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the cyclo(RGDyC) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

cyclo(RGDyC) is a cyclic pentapeptide that contains the αVβ3 integrin-binding sequence arginine-glycine-aspartate (RGD).^{1,2} It has been used to generate doxorubicin-containing liposomes and fluorescent silicon nanoparticles to target tumors and ocular neovascularization, respectively.

References

1. Amin, A., Mansourian, M., Koning, G.A., *et al.* Development of a novel cyclic RGD peptide for multiple targeting approaches of liposomes to tumor region. *J. Control Release* **220(Pt A)**, 308-315 (2015).
2. Tang, M., Ji, X., Xu, H., *et al.* Photostable and biocompatible fluorescent silicon nanoparticles-based theranostic probes for simultaneous imaging and treatment of ocular neovascularization. *Anal. Chem.* **90(13)**, 8188-8195 (2018).

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