

PRODUCT INFORMATION



Cyclo(RGDfC) (trifluoroacetate salt)

Item No. 35172

Formal Name: cyclo(L-arginylglycyl-L- α -aspartyl-D-phenylalanyl-L-cysteinyl), trifluoroacetate salt

Synonyms: c(RGDfC), cyclo(Arg-Gly-Asp-D-Phe-Cys)

MF: C₂₄H₃₄N₈O₇S • XCF₃COOH

FW: 578.6

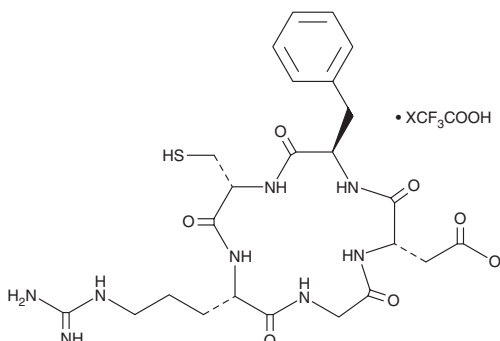
Purity: $\geq 90\%$

UV/Vis.: λ_{\max} : 204 nm

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(RGDfC) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the cyclo(RGDfC) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Cyclo(RGDfC) (trifluoroacetate salt) is soluble in ethanol and water.

Description

Cyclo(RGDfC) is a cyclic pentapeptide that contains the $\alpha\text{V}\beta 3$ integrin-binding sequence arginine-glycine-aspartate (RGD).^{1,2} It has been used in the synthesis of unimolecular micelles and ligand-decorated nanogels for cancer cell-targeted drug delivery.

References

1. Ryu, J.-H., Bickerton, S., Zhuang, J., *et al.* Ligand-decorated nanogels: Fast one-pot synthesis and cellular targeting. *Biomacromolecules* **13**(5), 1515-1522 (2012).
2. Xiao, Y., Hong, H., Javadi, A., *et al.* Multifunctional unimolecular micelles for cancer-targeted drug delivery and positron emission tomography imaging. *Biomaterials* **33**(11), 3071-3082 (2012).

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, 10/19/2022

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