

PRODUCT INFORMATION



RGD-targeted Proapoptotic Peptide

Item No. 42136

CAS Registry No.: 2243207-09-0

Formal Name: L-alanyl-L-cysteinyl-L- α -aspartyl-L-cysteinyl-L-arginylglycyl-L- α -aspartyl-L-cysteinyl-L-phenylalanyl-L-cysteinylglycylglycyl-D-lysyl-D-leucyl-L-alanyl-D-lysyl-D-leucyl-D-alanyl-D-lysyl-D-lysyl-D-leucyl-D-alanyl-D-lysyl-D-leucyl-D-alanyl-D-lysine, cyclic (2 $\rightarrow$ 10),(4 $\rightarrow$ 8)-bis(disulfide)

Peptide Sequence: A-cyclo[CD-cyclo[CRGDC]FC]GGKlAklAklaklak-OH

MF: C₁₁₆H₁₉₉N₃₅O₃₁S₄

FW: 2,708.3

Purity: $\geq$ 95%

Supplied as: A solid

Storage: -20°C

Stability: $\geq$ 4 years



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

Laboratory Procedures

RGD-targeted proapoptotic peptide is supplied as a solid. A stock solution may be made by dissolving the RGD-targeted proapoptotic peptide in the solvent of choice, which should be purged with an inert gas. RGD-targeted proapoptotic peptide is slightly soluble (0.1-1 mg/ml) in acetonitrile.

RGD-targeted proapoptotic peptide is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

RGD-targeted proapoptotic peptide is a peptide that contains an integrin-binding cyclic region (CDCRGDCFC) and a region similar to a pro-apoptotic heptadimer sequence (klaklaklaklak), both of which are found in similar peptides.¹

Reference

1. Gerlag, D.M., Borges, E., Tak, P.P., *et al.* Suppression of murine collagen-induced arthritis by targeted apoptosis of synovial neovasculature. *Arthritis Res.* **3**(6), 357-361 (2001).

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, 12/19/2024

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