

PRODUCT INFORMATION



DC-2 Peptide (trifluoroacetate salt)

Item No. 45287

Formal Name: (4S,7S,10S,13S,16S,19S,22S,25S,28S,31S,34S,37S,40S,43S)-1-amino-31-(2-amino-2-oxoethyl)-43-(((S)-1-(((S)-1-(((R)-4-((S)-1-(((S)-1-amino-3-(1H-indol-3-yl)-1-oxopropan-2-yl)amino)pyrrolidin-2-yl)-1-mercapto-3,4-dioxobutan-2-yl)amino)-3-(1H-imidazol-5-yl)-1-oxopropan-2-yl)amino)-3-(1H-imidazol-5-yl)-1-oxopropan-2-yl)carbonyl)-10,22,34,37-tetrakis(4-aminobutyl)-25-benzyl-7,13,19,28-tetrakis(3-guanidinopropyl)-4,16-diisobutyl-40-(2-(methylthio)ethyl)-2,5,8,11,14,17,20,23,26,29,32,35,38,41-tetradeca-oxo-3,6,9,12,15,18,21,24,27,30,33,36,39,42-tetradecaazapentatetracontan-45-oic acid, trifluoroacetate salt

H—Gly—Leu—Arg—Lys—Arg—Leu—Arg—Lys—Phe—Arg—
Asn—Lys—Lys—Met—Asp—His—His—Cys—Pro—Trp—NH₂
• XCF₃COOH

Peptide Sequence: H-GLRLRLRFRNLLMDHHCPW-NH₂

MF: C₁₁₅H₁₈₉N₄₃O₂₃S₂ • XCF₃COOH

FW: 2,606.2

Purity: ≥95%

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

DC-2 peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the DC-2 peptide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. DC-2 peptide (trifluoroacetate salt) is soluble (≥10 mg/ml) in ethanol and DMSO.

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 DC-2 peptide (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. DC-2 peptide (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

DC-2 peptide is a synthetic peptide that contains a DHHC palmitoylation motif.¹ It reduces H-Ras cell membrane localization in HeLa cells and increases ERK and Akt phosphorylation in SW480 colorectal cancer cells with high EGFR expression but not in SW620 colorectal cancer cells with low EGFR expression. DC-2 reduces viability of HeLa and MDA-MB-231 breast cancer cells at concentrations ranging from 20 to 100 μM.

Reference

1. Stillger, K., Platz-Baudin, E., Friedland, F., *et al.* First steps toward the design of peptides that influence the intracellular palmitoylation machinery. *ChemBiochem* **26**(10), e202500218 (2025).

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, 04/24/2026

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