

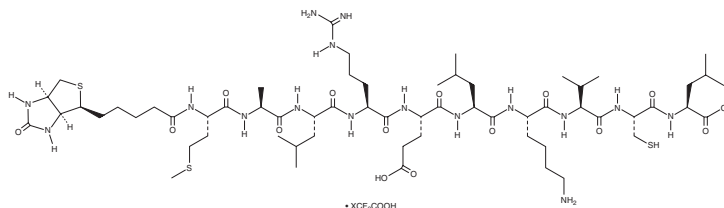
PRODUCT INFORMATION



Biotin-Rab-22A (1-10) (trifluoroacetate salt)

Item No. 44490

Formal Name: (2S,5R,8S,11S,14S,17S)-11-(4-aminobutyl)-17-((2S,5S,8S,11S)-2-(3-guanidinopropyl)-5-isobutyl-8-methyl-11-(2-(methylthio)ethyl)-4,7,10,13-tetraoxo-17-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)-3,6,9,12-tetraazaheptadecanamido)-2,14-diisobutyl-8-isopropyl-5-(mercaptomethyl)-4,7,10,13,16-pentaoxo-3,6,9,12,15-pentaazaicosanedioic acid, trifluoroacetate salt



Peptide Sequence: Biotin-MALRELKVCL-OH
MF: C₆₃H₁₀₉F₃N₁₆O₁₇S₃ • XCF₃COOH
FW: 1,515.8
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

Biotin-Rab-22A (1-10) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the biotin-Rab-22A (1-10) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Biotin-Rab-22A (1-10) (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO.

Description

Biotin-Rab-22A (1-10) is a biotinylated version of the GTPase Ras-related protein Rab-22A that corresponds to residues 1-10 of human and mouse Rab-22A.¹ It inhibits RhoA-GTP increases induced by Rab-22A (1-38) or an oncogenic Rab-22A fusion protein Rab-22A-NeoF1 in 143B whole cell lysates.

Reference

1. Liao, D., Zhong, L., Yin, J., *et al.* Chromosomal translocation-derived aberrant Rab22a drives metastasis of osteosarcoma. *Nat. Cell Biol.* **22(7)**, 868-881 (2020).

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, 12/09/2025

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