

PRODUCT INFORMATION



Fyn Peptide (410-430) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt)
Item No. 37507

Formal Name: (3S,6S)-N¹-((3S,6S,9S,15S,18S,21S,24S,27S,30S,35S)-37-amino-15-(3-amino-3-oxopropyl)-6-(4-aminobutyl)-3-benzyl-18-(3-((diaminomethylene)amino)propyl)-1-((S)-1-(hydroxy-L-isoleucyl)pyrrolidin-2-yl)-30-(hydroxyamino)-27-(4-hydroxybenzyl)-24-((R)-1-hydroxyethyl)-9,21-dimethyl-1,4,7,10,13,16,17,22,25,28,29,33,34,37-tetradecaoxo-2,5,8,11,14,19,20,23,26-nonaazaheptatriacontan-35-yl)-N⁹-((5S,8S,11S,14S,17S,18S)-11-(3-((diaminomethylene)amino)propyl)-5,14-diisobutyl-8,18-dimethyl-1,4,7,12,13,16-hexaoxo-3,6,9,10,15-pentaazaicosan-17-yl)-3,6-bis(hydroxyamino)-4,5-dioxononanediamide, trifluoroacetate salt

H—Gly—Leu—Ala—Arg—Leu—Ile—Glu—Asp—Asn—Glu—
Tyr—Thr—Ala—Arg—Gln—Gly—Ala—Lys—Phe—Pro—
Ile—OH
• XCF₃COOH

Synonyms: Src (409-429) (human, rat), Src (408-428) (mouse)

Peptide Sequence: GLARLIEDNEYTARQGAKFPI-OH

MF: C₁₀₅H₁₆₈N₃₀O₃₂ • XCF₃COOH

FW: 2,362.7

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

Fyn peptide (410-430) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Fyn peptide (410-430) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Fyn peptide (410-430) is a peptide fragment of the Src family non-receptor tyrosine kinase Fyn containing Tyr420, which undergoes autophosphorylation upon activation of full-length Fyn.¹

Reference

1. Weir, M.E., Mann, J.E., Corwin, T., *et al.* Novel autophosphorylation sites of Src family kinases regulate kinase activity and SH2 domain-binding capacity. *FEBS Lett.* **590(8)**, 1042-1052 (2016).

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, 11/08/2023

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