

PRODUCT INFORMATION



S1H (trifluoroacetate salt)

Item No. 37515

Synonym:	Site 1-binding Helix	
Peptide Sequence:	Ac-YIPKEQKYSFLQNPQT-NH ₂	Ac-Tyr-Ile-Pro-Lys-Glu-Gln-Lys-Tyr-Ser-Phe-
MF:	C ₉₄ H ₁₄₁ N ₂₃ O ₂₇ • XCF ₃ COOH	
FW:	2,025.3	Leu-Gln-Asn-Pro-Gln-Thr-NH ₂
Purity:	≥95%	• XCF ₃ COOH
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

S1H (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the S1H (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

S1H is a synthetic peptide mimetic of human growth hormone amino acids 36-51 and an antagonist of the growth hormone receptor.¹ It adopts an α -helical structure in solution and decreases STAT5 phosphorylation induced by human growth hormone in SK-MEL-28 and MALME-3M melanoma cell lines, as well as by prolactin in IM-9 B cell lymphocytes, when used at concentrations of 100 and 200 nM.

Reference

1. Basu, R., Nahar, K., Kulkarni, P., *et al.* A novel peptide antagonist of the human growth hormone receptor. *J. Biol. Chem.* **296**, 100588 (2021).

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, 10/04/2023

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