

PRODUCT INFORMATION



SPC25 Peptide (human, bovine) (trifluoroacetate salt)

Item No. 44404

Formal Name: L-asparaginy-L-leucyl-L-leucyl-L-lysyl-L-leucyl-L-isoleucyl-L-alanyl-L- α -glutamyl-L-valine, trifluoroacetate salt
Synonyms: Asn-Leu-Leu-Lys-Leu-Ile-Ala-Glu-Val-OH, Spindle Pole Body Component 25
Peptide Sequence: NLLKLIAEV-OH H-Asn-Leu-Leu-Lys-Leu-Ile-Ala-Glu-Val-OH
MF: $C_{47}H_{85}N_{11}O_{13} \cdot XCF_3COOH$ $\cdot XCF_3COOH$
FW: 1,012.3
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years

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

Laboratory Procedures

Spindle pole body component 25 (SPC25) peptide (human, bovine) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the SPC25 peptide (human, bovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SPC25 Peptide (human, bovine) (trifluoroacetate salt) is soluble (≥ 10 mg/ml) in DMSO.

Description

SPC25 peptide is a peptide fragment of SPC25, which is a protein member of the NDC80 kinetochore complex and an effector of kinetochore-microtubule attachment during cell division.

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, 10/14/2025

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