

# PRODUCT INFORMATION



## IκBα (20-39) (human) (trifluoroacetate salt)

Item No. 44044

<b>Synonyms:</b>	NFKBIA, NF-κB Inhibitor α
<b>Peptide Sequence:</b>	LKKERLLDDRHDHSGLDMSMKD-OH
<b>MF:</b>	$C_{98}H_{167}N_{31}O_{35}S \cdot XCF_3COOH$
<b>FW:</b>	2,371.7
<b>Purity:</b>	≥95%
<b>Supplied as:</b>	A solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥4 years

H—Leu—Lys—Lys—Glu—Arg—Leu—Leu—Asp—Asp—Arg—

His—Asp—Ser—Gly—Leu—Asp—Ser—Met—Lys—Asp—OH

• XCF<sub>3</sub>COOH

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

### Laboratory Procedures

IκBα (20-39) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the IκBα (20-39) (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. IκBα (20-39) (human) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

IκBα (20-39) (human) (trifluoroacetate salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

IκBα (20-39) is a peptide fragment of the N-terminal signal-receiving domain of IκBα (Item No. 43691), which is an inhibitor of NF-κB and a member of the IκB family of proteins.<sup>1,2</sup> It contains several post-translational modification sites that regulate IκBα activity, including the phosphorylation sites Ser32 and Ser36, the ubiquitination sites Lys21 and Lys22, and the SUMOylation sites Lys21 and Lys38.<sup>2</sup>

### References

1. Majdalawieh, A. and Ro, H.-S. Regulation of IκBα function and NF-κB signaling: AEBP1 is a novel proinflammatory mediator in macrophages. *Mediators Inflamm.* 823821 (2010).
2. Wang, X., Peng, H., Huang, Y., *et al.* Post-translational modifications of IκBα: The state of the art. *Front. Cell Dev. Biol.* 8, 574706 (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, 10/02/2025

#### 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