

PRODUCT INFORMATION



Bak BH3 (72-87) (human) (trifluoroacetate salt)

Item No. 36909

Formal Name: glycyl-L-glutaminy-L-valylglycyl-L-arginyl-L-glutaminy-L-leucyl-L-alanyl-L-isoleucyl-L-isoleucylglycyl-L- α -aspartyl-L- α -aspartyl-L-isoleucyl-L-asparaginy-L-arginine, trifluoroacetate salt

Peptide Sequence: GQVGRQLAIIGDDINR

MF: C₇₂H₁₂₅N₂₅O₂₄ • XCF₃COOH

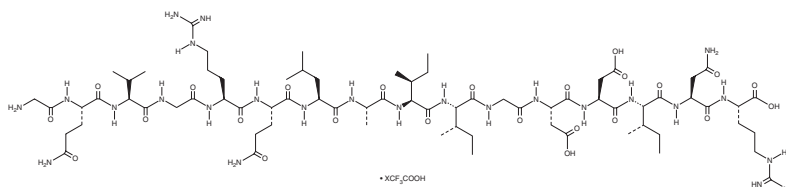
FW: 1,724.9

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

Bak BH3 (72-87) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Bak BH3 (72-87) (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Bak BH3 (72-87) (human) (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Bak BH3 (72-87) (human) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of Bak BH3 (72-87) (human) (trifluoroacetate salt) in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Bak BH3 (72-87) is a peptide that corresponds to amino acids 72-87 of the pro-apoptotic protein Bak.¹ It selectively binds to Bcl-xL over Bcl-2 (EC₅₀s = 174 and 3,020 nM, respectively). Bak BH3 (72-87) induces apoptosis in LLC-PK1 porcine kidney epithelial cells.²

References

1. Frey, V., Viaud, J., Subra, G., *et al.* Structure-activity relationships of Bak derived peptides: Affinity and specificity modulations by amino acid replacement. *Eur. J. Med. Chem.* **43**(5), 966-972 (2008).
2. Peherstorfer, E., Mayer, B., Boehm, S., *et al.* Effects of microinjection of synthetic Bcl-2 domain peptides on apoptosis of renal tubular epithelial cells. *Am. J. Physiol. Renal. Physiol.* **283**(1), F190-F196 (2002).

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, 12/21/2022

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