

PRODUCT INFORMATION

BimS BH3 (51-76) (human) (trifluoroacetate salt)

Item No. 36963

Formal Name:	L- α -aspartyl-L-methionyl-L-arginyl-L-prolyl-L- α -glutamyl-L-isoleucyl-L-tryptophyl-L-isoleucyl-L-alanyl-L-glutamyl-L- α -glutamyl-L-leucyl-L-arginyl-L-arginyl-L-isoleucylglycyl-L- α -aspartyl-L- α -glutamyl-L-phenylalanyl-L-asparaginyll-L-alanyl-L-tyrosyl-L-tyrosyl-L-alanyl-L-arginyl-L-arginine, trifluoroacetate salt	H—Asp—Met—Arg—Pro—Glu—Ile—Trp—Ile—Ala—Gln— Glu—Leu—Arg—Arg—Ile—Gly—Asp—Glu—Phe—Asn— Ala—Tyr—Tyr—Ala—Arg—Arg—OH
Synonyms:	DMRPEIWIAQELRRIGDEFNAYYARR-OH, Bim _s (51-76)	• XCF ₃ COOH
MF:	C ₁₄₅ H ₂₂₂ N ₄₄ O ₄₁ S • XCF ₃ COOH	
FW:	3,269.7	
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

BimS BH3 (51-76) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the BimS BH3 (51-76) (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. BimS BH3 (51-76) (human) (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BimS BH3 (51-76) (human) (trifluoroacetate salt) in these solvents is approximately 1, 2, and 3 mg/ml, respectively.

Description

BimS BH3 (51-76) is a peptide that corresponds to amino acids 51-76 of the short splice variant of the pro-apoptotic protein Bim (BimS), which contains the Bcl-2 homology 3 (BH3) motif.¹ It binds to myeloid cell leukemia-1 (Mcl-1), Bcl-xL, Bcl-2, and Bcl-W (IC₅₀s = 18, 10, 7, and 18 nM, respectively). BimS BH3 (51-76) induces cytochrome c release from mitochondria derived from *bcl-x*^{-/-} mouse embryonic fibroblasts (MEFs).

Reference

1. Lee, E.F., Fedorova, A., Zobel, K., *et al.* Novel Bcl-2 homology-3 domain-like sequences identified from screening randomized peptide libraries for inhibitors of the pro-survival Bcl-2 proteins. *J. Biol. Chem.* **284**(45), 31315-31326 (2009).

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