

Product Data Sheet

Product Name: Bim BH3, Peptide III

Catalog Number: AS-62278 (0.5 mg) Lot Number: See label on vial

Sequence: 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-OH (3-letter code) DMRPEIWIAQELRRIGDEFNAYYAR (1-letter code)

Molecular Weight: 3114.5

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 μl to 1 mg Bim BH3 peptide.

Storage: Bim BH3 peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: This Bim peptide belongs to the pro-apoptotic Bcl-2 family of proteins. Ref: Aoyagi, M. et al. *Prot. Sci.* **16**, 118 (2007).

Related Products:

NameCat #SizeBim BH3, Peptide IVAS-622791 mg(DMRPEIWIAQELRRIGDEFNAYYARR)

For Research Use Only