



Product Data Sheet

Product Name:	Noxa BH3, Peptide 1	
Catalog Number:	AS-62282 (1 mg)	Lot Number: See label on vial
Sequence:	H-Pro-Ala-Glu-Leu-Glu-Val-Glu-Cys-Ala-Thr-Gln-Leu-Arg-Arg-Phe-Gly-Asp-Lys-Leu-Asn-Phe-Arg-Gln-Lys-Leu-Leu-OH (3-letter code) PAELEVECATQLRRFGDKLNFRQKLL (1-letter code)	
Molecular Weight:	3075.6	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Noxa BH3 peptide is freely soluble in water.

Storage: Noxa 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 peptide belongs to the Bcl-2 family of proteins. Noxa gene encodes a Bcl-2 homology 3 (BH3)-only member of this family; it contains the BH3 region, not other BH domains. When ectopically expressed, Noxa undergoes BH3 motif-dependent localization to mitochondria, it interacts with anti-apoptotic Bcl-2 family members resulting in the activation of caspase-9. Ref: Oda, E. et al. *Science* **288**, 1053 (2000).

Related Products:

Name	Cat #	Size
Noxa A BH3 peptide, cell permeable rrrrrrrGAELPPEFAAQLRKIGDKVYC	AS-64087	1 mg

For Research Use Only