

## **Product Data Sheet**

Product Name: Proapoptotic Peptide, (klaklak)<sub>2</sub>

Catalog Number: AS-62199 (0.5 mg) Lot Number: See label on vial

Sequence: H-D-Lys-D-Leu-D-Ala-D-Lys-D-Leu-D-Ala-D-Lys-D-Leu-D-Ala-D-

Lys-D-Leu-D-Ala-D-Lys-NH2 (3-letter code)

klaklaklaklak-NH2 (1-letter code)

Molecular Weight: 1523.0

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Proapoptotic peptide is freely soluble in water.

Storage: Proapoptotic 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.

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

The synthetic peptides CGSPGWVRC, CGSPGWVRC-GG-D(KLAKLAK)2, and D(KLAKLAK)2 were obtained commercially (AnaSpec; San Jose, CA) to our specifications. In some binding experiments, as indicated, a synthetic cyclic peptide (sequence CARAC) served as an unrelated negative control. In Vitro Cell Viability and Cell Apoptosis Assays—Cells ( $2 \times 10^4$  cells/well) were seeded in 96-well plates for 24 h in 10% FBS in DMEM, incubated with increasing concentrations of the peptides CGSPGWVRC-GG-D(KLAKLAK)2 or CGSPGWVRC and D(KLAKLAK)2 in 80  $\mu$ l of 1% FBS in DMEM for 6 hours at 37 °C. After 6 h, cell viability was measured with a cell proliferation detection reagent according to the manufacturer's instructions (WST-1; Roche Applied Science)- Giordano, R. et al. J Biol Chem 283, 29447 (2008).

Soluble CGRRAGGSC-GG-D(KLAKLAK)<sub>2</sub>, CGRRAGGSC, and D(KLAKLAK)<sub>2</sub> peptides, and the unrelated control peptide CKGGRAKDC-GG-D(KLAKLAK)<sub>2</sub>, were synthesized to our specifications by AnaSpec (San Jose, CA). NCaP, MDA-PCa-2b cells (each at 3 x 10<sup>4</sup>/well), and EF43.*fgf*-4 cells (7) at 2 x 10<sup>4</sup>/well were seeded in triplicates and incubated in 96-well plates (Becton Dickinson, Franklin Lakes, NJ) for 24–72 h at 37°C, with serially increasing concentrations (10–100 μM) of CGRRAGGSC-GG-D(KLAKLAK)<sub>2</sub> peptide, CGRRAGGSC peptide alone, D(KLAKLAK)<sub>2</sub> peptide alone, or an equimolar mixture of the unconjugated peptides CGRRAGGSC and D(KLAKLAK)<sub>2</sub>. LNCaP cells were also exposed in parallel to increasing concentrations (20–100 μM) of CGRRAGGSC-GG-D(KLAKLAK)<sub>2</sub> and unrelated control peptides CKGGRAKDC-GG-D(KLAKLAK)<sub>2</sub> or CGSPGWVRC-GG-D(KLAKLAK)<sub>2</sub>, under the same conditions-Zurita, A. et al. *Cancer Res* 64, 435 (2004).

## **Published Citations:**

Zurita, A. et al. *Cancer Res* **64**, 435 (2004). Giordano, R. et al. *J Biol Chem* **283**, 29447 (2008). Nie, J. et al. *Stem Cells* **26**, 2735 (2008).

## Related Products:

Name Cat # Size
Pro-apoptotic Peptide, klaklakklaklak, 5-FAM-labeled AS-62206 1 mg
5-FAM-klaklakklaklak-NH2

For Research Use Only