



## Product Data Sheet

---

Product Name:	PCC (88 - 104)	
Catalog Number:	AS-60432-1 (1mg)	Lot Number: See label on vial
Sequence:	H-Lys-Ala-Glu-Arg-Ala-Asp-Leu-Ile-Ala-Tyr-Leu-Lys-Gln-Ala-Thr-Ala-Lys-OH (3-letter code) KAERADLIAYLKQATAK (1-letter code)	
Molecular Weight:	1890.2	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

**Peptide Reconstitution:** PCC (88-104) is freely soluble in water.

**Storage:** PCC (88-104) is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** PCC (88-104) is a 17-mer peptide fragment of pigeon cytochrome c that stimulates proliferative T cell responses. Ref: Wilson, DB. et al. *J. Immunol.* **163**, 6424 (1999).

**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.

Naive  $\text{CD4}^{+}$  T cells ( $5 \times 10^5$ ) were cultured for 24 h in 1 ml of RPMI in a 24-well plate with  $1.25 \times 10^5$  mitomycin C-treated P13.9 fibroblast cells stably expressing I-E<sup>k</sup>, CD80, and CD54, and preloaded with 0.001–1  $\mu\text{M}$  pPCC. Ref: Tarasenko, T. et al. *PNAS.* **104**, 11382 (2007).

### Published Citations:

Tarasenko, T. et al. *PNAS.* **104**, 11382 (2007).  
Stephan, TL. et al. *J. Immunol.* **183**. 5554 (2009).

*For Research Use Only*