

## **Product Data Sheet**

Product Name:	Phytochelatin 3, PC3	
Catalog Number:	AS-60790 (1 mg)	Lot Number: See label on vial
Sequence:	H- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys-Gly-OH (3-letter code) ( $\gamma$ E-C) <sub>3</sub> -G (1-letter code)	
Molecular Weight:	772.9	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Phytochelatin 3 peptide is freely soluble in water.

Storage: Phytochelatin 3 peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}$ C or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}$ C or lower.

Description: A glutathione-derived heavy metal-detoxifying peptide of higher plants consisting of 3 units of γGlu-Cys. Ref: Grill, E. et al. *Science* **230**, 674 (1985); Rauser, WE. *Plant Physiol.* **109**, 1141 (1995).

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.

Custom ordered phytochelatins (PC<sub>2</sub>, PC<sub>3</sub>, PC<sub>4</sub> and PC<sub>5</sub>) [PC<sub>n</sub>, ( $\gamma$ -Glu-Cys)<sub>n</sub>-Gly, where n = 2-5] were obtained from Anaspec, San Jose, CA, USA.Ten microliter aliquots of 8 mM stock solution of each standard (Cys, GSH,  $\gamma$ -EC, NAC, PC<sub>2</sub>, PC<sub>3</sub>, PC<sub>4</sub>, and PC<sub>5</sub>) were prepared using deionized water and stored in the dark at -20 °C. With the exception of NAC, appropriate portions of each stock were mixed together and further diluted with extraction buffer (6.3 mM DTPA with 0.1%, v/v, TFA) to create a series of seven working standards (S1–S7) with concentrations of Cys, GSH and  $\gamma$ -EC at 1, 2, 3, 4, 5, 7.5 and 10 pmol  $\mu$ L<sup>-1</sup> injected and PC<sub>2</sub>–PC<sub>5</sub> at 0.2, 0.4, 0.6, 0.8, 1.0, 1.5, and 2.0 pmol  $\mu$ L<sup>-1</sup> injected-Minocha, R. et al. *J. Chromatogr. A* **1207**, 72 (2008).

For quantification of PC<sub>n</sub>, respective standards ranging from PC<sub>1</sub> to PC<sub>5</sub> were obtained from Anaspec Inc. (Anaspec Inc., San Jose, CA, USA). All standard solutions were prepared and diluted in 1:1 acetonitrile:water solvent mixture. Separate stock solutions of 100 mg mL<sup>-1</sup> of each phytochelatin were prepared and stored at -80 °C. Aliquots of these solutions were mixed to obtain a 10 mg mL<sup>-1</sup> mixed working standard stock solution that was stored at -20 °C. Six-point calibration curves of mixed PC<sub>n</sub> analytes were prepared daily at 1, 10, 100, 250, 500, and 1000 µg mL<sup>-1</sup> concentrations using the 10 mg mL<sup>-1</sup> stock solution. The final volume was bought up to 0.5 mL using 1:1 acetonitrile:water solvent mixture and stored at -4 °C. Calibration curves were used for quantifying phytochelatins in the experimental plant samples-<u>Andra, SS. et al.</u> <u>Environ. Pollut. **157**, 2173 (2009).</u>

## **Published Citations:**

Kang, SH. et al. *Angewandte Chemie* **47**, 5186 (2008). Mendoza-Cozatl, DG. et al. *Plant J.* **54**, 249 (2008). Miao, AJ. & WX. Wang *Environ. Sci. Technol.* **41**, 1777 (2007). Minocha, R. et al. *J. Chromatogr. A* **1207**, 72 (2008). Zeng, X. et al. *Environ. Exp. Botany* **66**, 242 (2008). Andra, SS. et al. *Environ. Pollut.* **157**, 2173 (2009).

## **Related Products:**

Name Phytochelatin 2, PC2 (yE-C) <sub>2</sub> -G	<b>Cat #</b> AS-60791	<b>Size</b> 1 mg
Phytochelatin 4, PC4 (yE-C) <sub>4</sub> -G	AS-60789	1 mg
Phytochelatin 5, PC5 (yE-C) <sub>5</sub> -G	AS-61190	1 mg
Phytochelatin 6, PC6 (yE-C) <sub>6</sub> -G	AS-61191	1 mg

For Research Use Only