



service@anaspec.com • www.anaspec.com

# Phytochelatins

ISO 9001:2000 certified • 2149 O'Toole Ave. • San Jose, CA 95131 • 800-452-5530 • Fax 408-452-5059

## Phytochelatins

Product	MW	Size	Catalog#	US\$
<b>Phytochelatin, PC2</b> ( $\gamma$ E-C) <sub>2</sub> -G H- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys-Gly-OH A glutathione-derived heavy metal-detoxifying peptide of higher plants consisting of 2 units of $\gamma$ Glu-Cys. Ref: Grill, E. et al. <i>Science</i> <b>230</b> , 674 (1985); Rauser, WE. <i>Plant Physiol.</i> <b>109</b> , 1141 (1995).	540.6	1 mg	60791	\$90
<b>Phytochelatin, PC3</b> ( $\gamma$ E-C) <sub>3</sub> -G H- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys-Gly-OH A glutathione-derived heavy metal-detoxifying peptide of higher plants consisting of 3 units of $\gamma$ Glu-Cys. Ref: Grill, E. et al. <i>Science</i> <b>230</b> , 674 (1985); Rauser, WE. <i>Plant Physiol.</i> <b>109</b> , 1141 (1995).	772.9	1 mg	60790	\$95
<b>Phytochelatin, PC4</b> ( $\gamma$ E-C) <sub>4</sub> -G H- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys-Gly-OH A glutathione-derived heavy metal-detoxifying peptide of higher plants consisting of 4 units of $\gamma$ Glu-Cys. Ref: Grill, E. et al. <i>Science</i> <b>230</b> , 674 (1985); Rauser, WE. <i>Plant Physiol.</i> <b>109</b> , 1141 (1995).	1004.1	1 mg	60789	\$110
<b>Phytochelatin, PC5</b> ( $\gamma$ E-C) <sub>5</sub> -G H- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys-Gly-OH A glutathione-derived heavy metal-detoxifying peptide of higher plants consisting of 5 units of $\gamma$ Glu-Cys.	1236.4	1 mg	61190	\$120
<b>Phytochelatin, PC6</b> ( $\gamma$ E-C) <sub>6</sub> -G H- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys- $\gamma$ -Glu-Cys-Gly-OH A glutathione-derived heavy metal-detoxifying peptide of higher plants consisting of 6 units of $\gamma$ Glu-Cys.	1468.6	1 mg	61191	\$130