

## **Product Data Sheet**

Product Name:	Sauvagine	
Catalog Number:	AS-24475 (0.5 mg)	Lot Number: See label on vial
Sequence:	Pyr-Gly-Pro-Pro-IIe-Ser-IIe-Asp-Leu-Ser-Leu-Glu-Leu-Leu-Arg-Lys-Met- IIe Glu-IIe-Glu-Lys-Gln-Glu-Lys-Glu-Lys-Gln-Gln-Ala-Ala-Asn-Asn-Arg- Leu-Leu-Leu-Asp-Thr-IIe-NH2 (3-letter code) Pyr-GPPISIDLSLELLRKMIEIEKQEKEKQQAANNRLLLDTI-NH2 (1-letter code)	
Molecular Weight:	4599.4	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Initially, try to dissolve a small amount of Sauvagine peptide in water. If peptide does not dissolve in water, then use 1% NH<sub>4</sub>OH.

Storage: Sauvagine 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: A hypotensive and diuretic peptide. Originally isolated from the skin of the frog, Phyllomedusa sauvagei. It affects diuresis in the cardiovascular system and causes the release of ACTH and endorphins. Ref: Montecucchi, PC. et al. Int. *J. Pept. Protein Res.* **20**, 139 (1982); Erspamer, V. et al. *Trends Pharmacol. Sci.* **20**, 391 (1980).

## Published Citations:

Zheng, J. et al. Anal. Chem. 78, 1535 (2006).

Related Products:

Name Deltorphin A YmFHLMD-NH2 Cat # Size AS-62682 1 mg

For Research Use Only