



Product Data Sheet

Product Name:	Melanin Concentrating Hormone, human, mouse, rat	
Catalog Number:	22907 (0.5mg) 22908 (1mg)	Lot Number: See label on vial
Sequence:	H-Asp-Phe-Asp-Met-Leu-Arg-Cys-Met-Leu-Gly-Arg-Val-Tyr-Arg-Pro-Cys-Trp-Gln-Val-OH(Disulfidebridge:7-16) (3-letter code) DFDMLRCMLGRVYRPCWQV (Disulfide bridge: 7-16) (1-letter code)	
Molecular Weight:	2386.9	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Initially try to dissolve a small amount of Melanin Concentrating Hormone in water. If the peptide does not dissolve, use 1% NH_4OH .

Storage: Melanin Concentrating Hormone 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.

Published Citations:

Presse, F. et al. *Neurosci.* **71**, 735 (1996).
Vaughan, JM. et al. *Endocrinol.* **125**, 1660 (1989).

For Research Use Only