

Technical Data Sheet

Product Name	ACTH (1 - 17) SYSMEHFRWGKPVGKKR
Size	1 mg
Catalog #	AS-60708
Purity	% Peak Area By HPLC \geq 95%
Description	ACTH (1-17), and aMSH (a-Melanotropin), both derived from POMC (proopiomelanocortin) are involved in melanogenesis. However, ACTH (1-17) has been found to be more potent in melanogenesis in human melanocytes. Both ACTH (1-17) and aMSH also increase dendricity, and proliferation in follicular melanocytes.
Storage	-20°C
References	Kauser, S. et al. <i>FASEB J</i> 20 , 882 (2006) Tsatmali, M. et al. <i>Cell Mol Bio</i> 45 , 1029 (1999)
Molecular Weight	2093.4
Sequence (One-Letter Code)	SYSMEHFRWGKPVGKKR
Sequence (Three-Letter Code)	H - Ser - Tyr - Ser - Met - Glu - His - Phe - Arg - Trp - Gly - Lys - Pro - Val - Gly - Lys - Lys - Arg - OH