

Product Data Sheet

Product Name:	Kemptide		
Catalog Number:	AS-22594 (5 mg) AS-22595 (25 mg)	Lot Number: See label on vial	
Sequence:	H-Leu-Arg-Arg-Ala-Ser-Leu-Gly-OH (3-letter code) LRRASLG (1-letter code)		
Molecular Weight:	771.9		
Peptide Purity:	>95%		
Appearance:	Lyophilized white powder		

Peptide Reconstitution: Kemptide is freely soluble in water.

Storage: Kemptide 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: Kemptide is a phosphate acceptor peptide that serves as a synthetic substrate for PKA (Km = 16 μ M). The corresponding fluorescent and biotinylated peptides are also proven to be good substrates for PKA. Ref: Pfister, TD. and KB. Storey, *Insect Biochem. Mol. Biol.* **32**, 505 (2002); Zelada, A. et al. *Eur. J. Biochem.* **252**, 245 (1998).

Related Products:

Name	Cat #	Size
Kemptide [LRRASLG], 5-FAM labeled	AS-29933-1	1 mg
5-FAM-LRRASLG	AS-29933-5	5 mg

For Research Use Only