



Product Data Sheet

Product Name: ERKtide

Catalog Number: AS-60211-5 (5 mg) **Lot Number:** See label on vial

Sequence: H-Ala-Thr-Gly-Pro-Leu-Ser-Pro-Gly-Pro-Phe-Gly-Arg-Arg-OH
(3-letter code)
ATGPLSPGPFGRR (1-letter code)

Molecular Weight: 1313.5

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: ERKtide peptide is freely soluble in H₂O.

Storage: ERKtide 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: ERKtide is a peptide substrate for ERK2. Extracellular regulated protein kinase 2 (ERK2) is a eukaryotic protein kinase whose activity is regulated by mitogenic stimuli. Ref: Prowse, CN. et al. *Biochem.* **39**, 6258 (2000); Waas, W. et al. *Biochem.* **42**, 12273 (2003).

Related Products:

Name	Cat #	Size
Erktide (IPTTPITTTYFFFK)	AS-61777	1 mg

For Research Use Only