



## Product Data Sheet

---

**Product Name:** S6 Kinase Substrate (229-239)

**Catalog Number:** AS-27191 (1 mg)                      Lot Number: See label on vial  
AS-27192 (5 mg)

**Sequence:** H-Ala-Lys-Arg-Arg-Arg-Leu-Ser-Ser-Leu-Arg-Ala-OH (3-letter code)  
AKRRRLSSLRA (1-letter code)

**Molecular Weight:** 1313.6

**Peptide Purity:** >95%

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** S6 Kinase peptide is freely soluble in water.

**Storage:** S6 Kinase peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Reference:** House, C. et al. *J. Biol. Chem.* **262**, 772 (1987).

**Related Products:**

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
S6 Kinase Substrate (229-239), Amide, Biotinylated	AS-27193	1 mg
Biotin-AKRRRLSSLRA-NH <sub>2</sub>		

*For Research Use Only*