



## Product Data Sheet

---

<b>Product Name:</b>	CREBtide, 5-FAM labeled	
<b>Catalog Number:</b>	AS-60436-1 (1 mg) AS-60436-5 (5 mg)	Lot Number: See label on vial
<b>Sequence:</b>	5-FAM-Lys-Arg-Arg-Glu-Ile-Leu-Ser-Arg-Arg-Pro-Ser-Tyr-Arg-OH (3-letter code) 5-FAM-KRREILSRRPSYR (1-letter code)	
<b>Molecular Weight:</b>	2075.3	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized yellow powder	

**Peptide Reconstitution:** CREBtide is freely soluble in water.

**Storage:** CREBtide 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.

**Description:** CREBtide (60510-1) is a synthetic substrate for PKA ( $K_m = 3.9 \mu\text{M}$ ). This peptide is based on the phosphorylation sequence in  $\delta$ -CREB (cAMP response element binding protein).  
Ref: Bok, J. et al. *J. Neurosci.* **23**, 777 (2003); Colbran, J.L. et al. *Biochem. Cell Biol.* **70**, 1277 (1992). Yamagata, Y. et al. *J. Biol. Chem.* **266**, 15391 (1991).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
CREBtide, Phosphorylated, C-term K KRREILSRRPS-pY-RK	AS-27018	1 mg
CREBtide KRREILSRRPSYR	AS-60510-5	5 mg

*For Research Use Only*