# **CREBtide**

#### **PKA** substrate

This peptide, KRREILSRRPSYR, is an excellent substrate for many serine kinases, including PKA and PKC.

## **Ordering Information**

Order Online »

**BML-P195-0001** 1mg

Manuals, SDS & CofA

View Online »

### **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended. Store, as supplied,

at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

Long Term Storage -20°C

Shipping Blue Ice

### Regulatory Status RUO - Research Use Only

#### **Product Details**

**Appearance** White to off-white solid.

Formulation Lyophilized.

**Identity** Determined by MS.

**MW** 1718

Purity ≥95% (HPLC)

**Sequence** Lys-Arg-Glu-Ile-Leu-Ser-Arg-Pro-Ser-Tyr-Arg

**Solubility** Soluble in water.

**Source** Synthetic.

Technical Info / Product Notes Replacement for ADI-SPK-102.

Last modified: May 29, 2024

