



## Product Information Sheet

---

**Product Name:** Phosphopeptide Standard Kit  
**Size:** 6 Phosphopeptides at 10 µg each (6 vials)  
**Catalog Number:** AS-61145

<b>Phosphopeptides:</b>	<b>MW</b>
FQ-pS-EEQQQTEDELQDK (Bovine β-Casein, monophospho cat# AS-61146)	2062.0
KRP-pS-QRHGSKY-NH <sub>2</sub> (UOM9, Phosphorylated PKC Substrate-1, cat# AS-20294)	1422.5
TRDIYETDpYYRK (Kinase Domain of Insulin Receptor, cat# AS-20274)	1702.8
TRDIpYETDpYpYRK (Kinase Domain of Insulin Receptor, cat# AS-20272)	1862.8
DLDVPIPGRFDRRV-pS-VAAE (PKA Regulatory Subunit II Substrate)	2192.4
SFVLNPTNIGM-pS-KSSQGHVTK (DAM1 [221-241])	2312.6

**HPLC Purity:** > 95%

**Storage:** -20°C

### Instructions:

Dissolve each phosphopeptide in 50 µl of 50% Acetonitrile. Spot 2-3 µl onto a MALDI target plate.

### References:

- Posewitz, MC. and P. Tempst, *Anal. Chem.* **71**, 2883 (1999)
- Show, W. et al. *J. Am. Soc. Mass Spectrom.* **11**, 273 (2000)
- Kim, J. et al. *J. Mass Spectrom.* **39**, 208 (2004)
- Liu, S. et al. *Rapid Commun. Mass Spectrom* **19**, 2747 (2005).

