



## SensoLyte® Protein Kinase Substrate Sampler Kit *\*Biotinylated\**

<b>Catalog #</b>	<b>71200</b>
<b>Unit Size</b>	1 kit

This kit contains ten biotinylated protein kinase peptide substrates and two phosphorylated peptide controls. It is designed to facilitate the identification of the optimum peptide substrate for your protein kinase of interest.

### USA and Canada Ordering Information

#### **AnaSpec Corporate Headquarter**

2149 O'Toole Ave.  
San Jose, CA 95131  
Toll-Free: 800-452-5530  
Tel: 408-452-5055  
Fax: 408-452-5059  
E-mail: [service@anaspec.com](mailto:service@anaspec.com)  
Internet: [www.anaspec.com](http://www.anaspec.com)

### International Ordering Information

A list of international distributors is available at [www.anaspec.com](http://www.anaspec.com).

## **INTRODUCTION**

Protein kinases play critical roles in signal transduction and are important drug-screening targets. The identification of an appropriate kinase substrate has proven to be critical in designing kinase assays. The SensoLyte® Protein Kinase Substrate Sampler Kit provides ten biotinylated protein kinase substrates along with two phosphorylated peptide controls (refer to Table 1 for peptide sequences). These substrates can be utilized to design ELISA-based kinase assays, facilitating the identification of an appropriate peptide substrate for the protein kinase of interest. All the components in this kit are available individually and in bulk packaging.

## **KIT COMPONENTS, STORAGE AND HANDLING**

*Note: Store all peptides at -20 °C, and keep away from light.*

**Component A:** Protein kinase substrate, biotinylated, (50 µL 1 mM DMSO solution X 12 vials)

Table 1. Protein kinase peptide substrate sequences.

Substrate No.	Catalog No.	Kinase	Kinase Substrate Peptide Sequence	FW
SB 1	29945-1	AKT/PKB/Rac-protein kinases	<b>Biotin</b> -ARKRERTYSFGHHA <sup>1:2</sup>	1942.2
SB 2	29915-1	Glycogen Synthase (1-8)	<b>Biotin</b> -PLSRTLVAACK-NH <sub>2</sub> <sup>3-5</sup>	1495.9
SB 3	29930-1	Ca <sup>2+</sup> -calmodulin-dependent protein kinase	<b>Biotin</b> -KKALRRQETVDAL <sup>6-8</sup>	1754.1
SB 4	29975-1	Ca <sup>2+</sup> -calmodulin-dependent protein kinase II	<b>Biotin</b> -KKALHRQETVDAL <sup>9:10</sup>	1735.0
SB 5	29957-1	CaM kinase II	<b>Biotin</b> -MHRQETVDCLK-NH <sub>2</sub> <sup>11:12</sup>	1584.9
SB 6	29918-1	MARCKS Protein (154-165)	<b>Biotin</b> -KKRFSFKKSFKL <sup>13-15</sup>	1770.2
SB 7	60513-1	Glycogen synthase kinase-3	<b>Biotin</b> -GRPRTSSFAEG <sup>2:16;17</sup>	1390.5
SB 8	29927-1	PKC $\mu$	<b>Biotin</b> -AALVRQMSVAFFFK <sup>18-20</sup>	1841.3
SB 9	60027-1	CDK5	<b>Biotin</b> -PKTPKKAKKL <sup>21;22</sup>	1365.8
SB 10	29966-1	MAP kinase (ERK1/ERK2 or p44/p42 MAPK)	<b>Biotin</b> -KRELVEPLTPSGEAPNQALLR-NH <sub>2</sub> <sup>23;24</sup>	2544.0
Control 1 for SB 1	29946-1	AKT/PKB/Rac-protein kinases	ARKRERTYS(PO <sub>3</sub> )FGHHA	1795.9
Control 2 for SB 3	29929-1	Ca <sup>2+</sup> -calmodulin-dependent protein kinase	KKALRRQE(PO <sub>3</sub> )TVDAL	1607.8

## **References**

1. T. Obata et al., *J.Biol.Chem.* 275, 36108-36115 (2000).

2. T. O. Chan, S. E. Rittenhouse, P. N. Tschlis, *Annu.Rev.Biochem.* 68, 965-1014 (1999).
3. D. R. Alexander, J. D. Graves, S. C. Lucas, D. A. Cantrell, M. J. Crumpton, *Biochem.J.* 268, 303-308 (1990).
4. C. House, R. E. Wettenhall, B. E. Kemp, *J.Biol.Chem.* 262, 772-777 (1987).
5. R. B. Pearson, J. R. Woodgett, P. Cohen, B. E. Kemp, *J.Biol.Chem.* 260, 14471-14476 (1985).
6. H. Takano, H. Fukushi, Y. Morishima, Y. Shirasaki, *Brain Res.* 962, 41-47 (2003).
7. R. E. Blair, S. B. Churn, S. Sombati, J. K. Lou, R. J. DeLorenzo, *Brain Res.* 851, 54-65 (1999).
8. P. M. Jones and S. J. Persaud, *Am.J.Physiol* 274, E708-E715 (1998).
9. M. Fluck, M. N. Waxham, M. T. Hamilton, F. W. Booth, *J.Appl.Physiol* 88, 352-358 (2000).
10. E. Yang and H. Schulman, *J.Biol.Chem.* 274, 26199-26208 (1999).
11. K. A. Kumar and P. P. Babu, *Neurosci.Lett.* 336, 121-125 (2003).
12. Y. Yamagata, A. J. Czernik, P. Greengard, *J.Biol.Chem* 266, 15391-15397 (1991).
13. K. Irie et al., *Pharmacol.Ther.* 93, 271-281 (2002).
14. F. Uberall et al., *J.Biol.Chem.* 272, 4072-4078 (1997).
15. P. J. Blackshear, *J.Biol.Chem.* 268, 1501-1504 (1993).
16. F. Dai et al., *Biochem.Biophys.Res.Commun.* 293, 1191-1196 (2002).
17. R. Meier, D. R. Alessi, P. Cron, M. Andjelkovic, B. A. Hemmings, *J.Biol.Chem.* 272, 30491-30497 (1997).
18. H. J. Kim et al., *J.Biol.Chem.* 278, 319-326 (2003).
19. K. Nishikawa, A. Toker, F. J. Johannes, Z. Songyang, L. C. Cantley, *J.Biol.Chem.* 272, 952-960 (1997).
20. I. Brandlin, T. Eiseler, R. Salowsky, F. J. Johannes, *J.Biol.Chem* 277, 45451-45457 (2002).
21. J. Lew, *Biochemistry* 42, 849-856 (2003).
22. P. Sharma et al., *J.Biol.Chem* 274, 9600-9606 (1999).
23. F. Sommer et al., *Int.J.Impot.Res.* 14, 217-225 (2002).
24. J. K. Klarlund, N. Khalaf, L. Kozma, M. P. Czech, *J.Biol.Chem* 268, 7646-7649 (1993).