

## BRSK2

(BR serine/threonine-protein kinase 2, Serine/threonine-protein kinase 29  
Serine/threonine kinase SAD-A, PEN11B, STK29)

- **CAT#:** BRSK2
- **Enzyme:** Human BRSK2
- **Substrate:** Peptide substrate, [KKLNRTLSEFAEPG], 5  $\mu$ M  
ATP 1  $\mu$ M
- **Reaction:**  

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.62
Ro 31-8220	77.6
GW 5074	6,300
H-89	815

- **Scientific Information:** Q8IWQ3 (<http://www.expasy.org/uniprot/Q8IWQ3>)

### Reference Compound IC50 for BRSK2

