

## NEK11

(Serine/threonine-protein kinase Nek11, NimA-related protein kinase 11, Never in mitosis A-related kinase 11)

- **CAT#:** NEK11
- **Enzyme:** Human NEK11
- **Substrate:** MBP (Myelin Basic Protein), 20  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**  

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

Compounds	IC50 (nM)
Staurosporine	570
Ro 31-8220	17,800
GW5074	2,000
Rottlerin	6,600

- **Scientific Information:** Q8NG66 (<http://www.expasy.org/uniprot/Q8NG66>)

### Reference compound IC50 for NEK11

