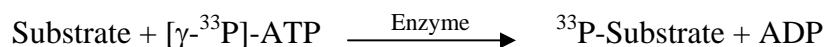


NEK7

(Serine/threonine-protein kinase Nek7, NimA-related protein kinase 7)

- **CAT#:** NEK7
- **Enzyme:** Human NEK7
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	19,500
PKR Inhibitor	6,500

- **Scientific Information:** Q8TDX7 (<http://www.expasy.org/uniprot/Q8TDX7>)

Reference compound IC50 for NEK7

