

TBK1

(Serine/threonine-protein kinase TBK1, TANK-binding kinase 1, T2K, NF-kappa-B-activating kinase, NAK)

- **CAT#:** TBK1
- **Enzyme:** Human TBK1
- **Substrate:** Peptide substrate, [KRRRAL[pS]VASLPGL], 20 μ M
ATP 10 μ 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	0.6
Ro 31-8220	530
GW 5074	5,800
Rottlerin	>20,000

- **Scientific Information:** Q9UHD2 (<http://www.expasy.org/uniprot/Q9UHD2>)

Reference compound IC50 for TBK1

