

ITK

(Tyrosine-protein kinase ITK/TSK, T-cell-specific kinase,
Tyrosine-protein kinase Lyk, Kinase EMT)

- **CAT#:** ITK
- **Enzyme:** Human ITK
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M
- **Reaction:**

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

Compounds	IC50 (nM)
Staurosporine	6.4
PP2	>20,000
AG1478	19,000
PDGFR I-III	>20,000
D-64406	3,000

- **Scientific Information:** Q08881 (<http://www.expasy.org/uniprot/Q08881>)

Reference compound IC50 for ITK

