

## FLT3

(FL cytokine receptor [Precursor], Tyrosine-protein kinase receptor FLT3, Stem cell tyrosine kinase 1, STK-1, Flk-2)

- **CAT#:** FLT3
- **Enzyme:** Human FLT3
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 5  $\mu$ 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	2.2
PP2	2,900
AG1478	320
PDGFR I-III	302
D-64406	56.5

- **Scientific Information:** P36888 (<http://www.expasy.org/uniprot/P36888>)

Reference compound IC50 for FLT3

