

## ABL1

(Proto-oncogene tyrosine-protein kinase ABL1, c-ABL, p150, Abelson murine leukemia viral oncogene homolog 1)

- **CAT#:** ABL1
- **Enzyme:** Human ABL1
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 20  $\mu$ M  
ATP 10  $\mu$ 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	60
PP2	140
AG1478	540
D-64406	670

- **Scientific Information:** P00519 (<http://www.expasy.org/uniprot/P00519>)

### Reference compound IC50 for ABL1

