

ABL1 (T315I)

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

- **CAT#:** ABL1 (T315I)
- **Enzyme:** Human ABL1 (T315I)
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 5 μ M
ATP 10 μ 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	24.2
AG1478	2,400
D-64406	2,600

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

Reference Compound IC50 for ABL1 (T315I)

