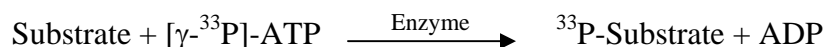


ABL2/ARG

(Tyrosine-protein kinase ABL2, Abelson murine leukemia viral oncogene homolog 2, Abelson-related gene protein, Tyrosine kinase ARG, ABLL)

- **CAT#:** ABL2
- **Enzyme:** Human ABL2
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	17
PP2	162
AG1478	580
D-64406	1,700

- **Scientific Information:** P42684 (<http://www.expasy.org/uniprot/P42684>)

Reference compound IC50 for ABL2/ARG

