

---

## ABL(ABL1)[T315I]

Product code 08-093

Full-length human ABL [2-1130(end) amino acids and T315I of accession number NP\_005148.2] was expressed as N-terminal His-tagged protein (126 kDa) using baculovirus expression system. His-tagged ABL[T315I] was purified by using Ni-NTA affinity chromatography.

Assay platform : Mobility Shift Assay

Substrate : ABLtide

ATP ( $\mu$ M) Km app / Bin : 4 / 5

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 6.4

IC50 at 1 mM ATP (nM) : 890

---