
ABL(ABL1)[E255K]

Product code: 08-094

Full-length human ABL [2-1130(end) amino acids and E255K of accession number NP_005148.2] was expressed as N-terminal His-tagged protein (126 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : ABLtide

ATP (μ M) Km app / Bin : 17 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 140

IC50 at 1 mM ATP (nM) : 4500
