
ABL(ABL1)

Product code: 08-001

Full-length human ABL [2-1130(end) amino acids 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 : 16 / 25
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 75
IC50 at 1 mM ATP (nM) : 1300
