
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

