
KIT[V654A]

Product code: 08-511

Human KIT, cytoplasmic domain [544-976(end) amino acids and V654A of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Srctide

ATP (μ M) Km app / Bin : 220 / 250

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 4.5

IC50 at 1 mM ATP (nM) : 8.2
