
PDGFR α (PDGFRA)

Product code 08-157

Human PDGFR α , cytoplasmic domain [550-1089(end) amino acids of accession number NP_006197.1] was expressed as N-terminal GST-fusion protein(89 kDa) using baculovirus expression system. GST-PDGFR α was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CSKtide

ATP (μ M) Km app / Bin : 28 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.30

IC50 at 1 mM ATP (nM) : 1.4
