
PDGFR β (PDGFRB)

Product code 08-158

Human PDGFR β , cytoplasmic domain [557-1106(end) amino acids of accession number NP_002600.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-PDGFR β was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CSKtide

ATP (μ M) K_{mapp} / B_{in} : 23 / 25

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 0.27

IC₅₀ at 1 mM ATP (nM) : 0.62
