
PDGFR α (PDGFRA) [D842V]

Product code 08-506

Human PDGFR α , cytoplasmic domain [550-1089(end) amino acids and D842V of accession number NP_006197.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-PDGFR α [D842V] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CSKtide

ATP (μ M) K_{mapp} / Bin : 21 / 25

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP Bin (nM) : 0.25

IC₅₀ at 1 mM ATP (nM) : 1.9
