
PDGFR α (PDGFRA)[V561D]

Product code: 08-507

Human PDGFR α , cytoplasmic domain [550-1089(end) amino acids and V561D of accession number NP_006197.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 35 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.32
IC50 at 1 mM ATP (nM)	: 1.6

