
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 |

