

---

---

## PDGFR $\alpha$ (PDGFRA)[D842V]

Product code: 08-506

Human PDGFR $\alpha$ , 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.

Assay platform	: MSA
Substrate	: CSKtide
ATP ( $\mu$ M) Km app / Bin	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.25
IC50 at 1 mM ATP (nM)	: 1.9

