
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.

Assay platform	: MSA
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 23 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.27
IC50 at 1 mM ATP (nM)	: 0.62

