
IGF1R

Product code: 08-141

Human IGF1R, cytoplasmic domain [959-1367(end) amino acids of accession number NP_000866.1] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : IRS1

ATP (μ M) Km app / Bin : 63 / 75

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 40

IC50 at 1 mM ATP (nM) : 150
