
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. GST-IGF1R was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : IRS1

ATP (μ M) K_{mapp} / B_{in} : 63 / 75

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 40

IC₅₀ at 1 mM ATP (nM) : 150
