
INSR

Product code 08-142

Human INSR, catalytic domain [1005-1310 amino acids of accession number NP_000199.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-INSR was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : IRS1

ATP (μ M) K_{mapp} / B_{in} : 58 / 50

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 12

IC₅₀ at 1 mM ATP (nM) : 70
