
IRR(INSRR)

Product code: 08-143

Human IRR, cytoplasmic domain [953-1297(end) amino acids of accession number NP_055030.1] was expressed as N-terminal GST-fusion protein (66 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : IRS1

ATP (μ M) Km app / Bin : 64 / 75

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 15

IC50 at 1 mM ATP (nM) : 98
