
DDR2

Product code: 08-114

Human DDR2, cytoplasmic domain [422-855(end) amino acids of accession number NP_006173.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : IRS1

ATP (μ M) Km app / Bin : 38 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.2

IC50 at 1 mM ATP (nM) : 0.77
