
DDR1

Product code: 08-113

Human DDR1, cytoplasmic domain [444-876(end) amino acids of accession number NP_001945.3] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : IRS1

ATP (μ M) Km app / Bin : 94 / 100

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 4.0

IC50 at 1 mM ATP (nM) : 3.1
