
RON(MST1R)

Product code: 08-152

Human RON, cytoplasmic domain [979-1400(end) amino acids of accession number NP_002438.2] was expressed as N-terminal GST-fusion protein (75kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Srctide

ATP (μ M) Km app / Bin : 27 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 37

IC50 at 1 mM ATP (nM) : 550
