
RON(MST1R)

Product code: 08-152

Human RON, cytoplasmic domain [979-1400(end) amino acids of accession number NP_002438.1] 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

