
ROS(ROS1)

Product code: 08-163

Human ROS, cytoplasmic domain [1883-2347(end) amino acids of accession number NP_002935.2] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : IRS1

ATP (μ M) Km app / Bin : 37 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 0.4
IC50 at 1 mM ATP (nM) : 1.0
