

---

## ***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. GST-ROS was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : IRS1

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 37 / 50  
Metal : Mg  
Reference compound : Staurosporine  
IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 0.4  
IC<sub>50</sub> at 1 mM ATP (nM) : 1.0

---