

---

---

# 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 ( $\mu$ 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

