

---

# YES(YES1)

Product code: 08-175

Full-length human YES [1-543(end) amino acids of accession number NP\_005424.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Src tide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 13 / 10

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 2.4

IC<sub>50</sub> at 1 mM ATP (nM) : 23

---