

---

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

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 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

---