

---

---

# 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                   | : Srctide       |
| ATP ( $\mu$ M) Km app / Bin | : 13 / 10       |
| Metal                       | : Mg            |
| Reference compound          | : Staurosporine |
| IC50 at ATP Bin (nM)        | : 2.4           |
| IC50 at 1 mM ATP (nM)       | : 23            |

