

---

# FER

Product code 08-139

Full-length human FER [1-822(end) amino acids of accession number NP\_005237.1] was expressed as N-terminal GST-fusion protein (122 kDa) using baculovirus expression system. GST-FER was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 26 / 25

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 1.0

IC<sub>50</sub> at 1 mM ATP (nM) : 12

---