

---

# FES

Product code: 08-140

Full-length human FES [1-413, 416-822(end) amino acids of accession number NP\_001996.1] was expressed as N-terminal GST-fusion protein (120 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 43 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.9
IC50 at 1 mM ATP (nM)	: 25

