

---

# Haspin(GSG2)

Product code 05-111

Full-length human Haspin [1-798(end) amino acids of accession number NP\_114171.2] was expressed as N-terminal GST-fusion protein (116 kDa) using baculovirus expression system. GST-Haspin was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : Histone H3 peptide

ATP ( $\mu$ M) Km app / Bin : 140 / 150  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 5.8  
IC50 at 1 mM ATP (nM) : n.a.

---