
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.

Assay platform	: MSA
Substrate	: Histone H3 peptide
ATP (μ 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.
