

---

# BUB1/BUB3

Product code: 05-187

Full-length human BUB1 [1-1085(end) amino acids of accession number NP\_004327] was co-expressed as N-terminal GST-fusion protein (149 kDa) with DYKDDDDK tagged BUB3 [1-328 (end) amino acids of accession number NP\_004716] using baculovirus expression system. GST-BUB1 was purified by using glutathione sepharose chromatography.

Assay platform : MSA  
Substrate : H2A peptide

ATP ( $\mu$ M) Km app / Bin : 2.9 / 5  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 300  
IC50 at 1 mM ATP (nM) : n.a.

---