
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 (μ 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.
