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

Substrate : H2A peptide

: Mobility Shift Assay

ATP (μ M) Kmapp / Bin : 2.9 / 5 Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 300

IC50 at 1 mM ATP (nM): n.a.