
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.

Assay platform : MSA
Substrate : H2A peptide

ATP (μ M) K_m app / B_{in} : 2.9 / 5
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 300
IC₅₀ at 1 mM ATP (nM) : n.a.
