
AurC(AURKC)

Product code 05-103

Full-length human AurC [1-275(end) amino acids of accession number NP_003151.2] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system. GST-AurC was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : Kemptide

ATP (μ M) Kmapp / Bin : 24 / 25
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 3.1
IC50 at 1 mM ATP (nM) : 18
