
GSK3β(GSK3B)

Product code 04-141

Full-length human GSK3β [1-420(end) amino acids of accession number NP_001139628.1] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-GSK3β was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : CREBtide-p

ATP (μM) Km app / Bin : 9.1 / 10
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 9.2
IC50 at 1 mM ATP (nM) : 240
