
GSK3α(GSK3A)

Product code 04-140

Full-length human GSK3α [1-483(end) amino acids of accession number NP_063937.2] was expressed as N-terminal GST-fusion protein (78 kDa) using baculovirus expression system. GST-GSK3α was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CREBtide-p

ATP (μM) Km app / Bin : 12 / 10

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 15

IC50 at 1 mM ATP (nM) : 400
