
GSK3 β (GSK3B)

Product code: 04-141

Full-length human GSK3 β [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.

Assay platform	: MSA
Substrate	: CREBtide-p
ATP (μ M) 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
