

---

---

## **GSK3 $\alpha$ (GSK3A)**

Product code: 04-140

Full-length human GSK3 $\alpha$  [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.

Assay platform	: MSA
Substrate	: CREBtide-p
ATP ( $\mu$ M) Km app / Bin	: 12 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 15
IC50 at 1 mM ATP (nM)	: 400

---

---