

---

## CGK2(PRKG2)

Product code 01-143

Full-length human CGK2 [1-762(end) amino acids of accession number NP\_006250.1] was expressed as N-terminal GST-fusion protein (114 kDa) using baculovirus expression system. GST-CGK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Kemptide

ATP ( $\mu$ M) Km app / Bin : 24 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.88

IC50 at 1 mM ATP (nM) : n.a.

---