

---

## **CDC7/ASK**

Product code 05-109

Full-length human CDC7 [1-574(end) amino acids of accession number NP\_003494.1] was co-expressed as N-terminal GST-fusion protein (92 kDa) with Dbf4(ASK) [1-674(end) amino acids of accession number NP\_006707.1] using baculovirus expression system. GST-CDC7 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : MCM2 peptide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 2.8 / 5

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 16

IC<sub>50</sub> at 1 mM ATP (nM) : 1600

---