

---

# DYRK1A

Product code 04-130

Full-length human DYRK1A [1-763(end) amino acids of accession number NP\_001387.2] was expressed as N-terminal GST-fusion protein (112 kDa) using baculovirus expression system. GST-DYRK1A was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : DYRKtide-F

ATP ( $\mu$ M) Km app / Bin : 16 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 7.8

IC50 at 1 mM ATP (nM) : 120

---