

---

# DYRK2

Product code 04-132

Full-length human DYRK2 [1-528(end) amino acids of accession number NP\_003574.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-DYRK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : DYRKtide-F

ATP ( $\mu$ M) Km app / Bin : 7.7 / 10

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 130

IC50 at 1 mM ATP (nM) : 6300

---