

---

# 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.

Assay platform	: MSA
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

