

---

---

# DYRK1B

Product code: 04-131

Full-length human DYRK1B [1-629(end) amino acids of accession number NP\_004705.1] was expressed as N-terminal GST-fusion protein (96 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DYRKtide-F
ATP ( $\mu$ M) Km app / Bin	: 59 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: 32

