

---

# DYRK3

Product code: 04-133

Full-length human DYRK3 [1-588(end) amino acids of accession number NP\_003573.2] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DYRKtide-F
ATP ( $\mu$ M) Km app / Bin	: 6.8 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 17
IC50 at 1 mM ATP (nM)	: 1200

