
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. GST-DYRK3 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform : Mobility Shift Assay

Substrate : DYRKtide-F

ATP (μ 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
