
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 (μ 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
