
DYRK1A

Product code: 04-130

Full-length human DYRK1A [1-763(end) amino acids of accession number NP_001387.2] was expressed as N-terminal GST-fusion protein (112 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DYRKtide-F
ATP (μ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 7.8
IC50 at 1 mM ATP (nM)	: 120

