
NuaK1

Product code: 02-126

Full-length human NuaK1 [1-661(end) amino acids of accession number NP_055655.1] was expressed as N-terminal GST-fusion protein (102 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : CHKtide

ATP (μ M) Km app / Bin : 59 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 1.0
IC50 at 1 mM ATP (nM) : 3.4
