
NuaK2

Product code: 02-127

Full-length human NuaK2 [1-628(end) amino acids of accession number NP_112214.3] was expressed as N-terminal GST-fusion protein (98kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : CHKtide

ATP (μ M) Km app / Bin : 26 / 25
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 2.2
IC50 at 1 mM ATP (nM) : 2.3
