
NuaK2

Product code 02-127

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

Assay platform : Mobility Shift Assay

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
