
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. GST-NuaK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

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
