
DAPK1

Product code 02-134

Human DAPK1, catalytic domain [1-289 amino acids of accession number NP_004929.1] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-DAPK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : DAPK1tide

ATP (μ M) K_{mapp} / B_{in} : 1.1 / 1

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 2.9

IC₅₀ at 1 mM ATP (nM) : 490
