
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.

Assay platform : MSA
Substrate : DAPK1tide

ATP (μ M) K_m app / 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
