
ACK(TNK2)

Product code: 08-196

Human ACK, catalytic domain [110-476 amino acids of accession number NP_005772.3] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : WASP peptide

ATP (μ M) Km app / Bin : 97 / 100
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 3.2
IC50 at 1 mM ATP (nM) : 3.8
