
LTK

Product code: 08-106

Human LTK, catalytic domain [498-796 amino acids of accession number NP_002335.2] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Src tide
ATP (μ M) K_m app / B_{in}	: 49 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC ₅₀ at ATP B_{in} (nM)	: 2.0
IC ₅₀ at 1 mM ATP (nM)	: 7.1