

---

---

# 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	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 49 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.0
IC50 at 1 mM ATP (nM)	: 7.1

