

---

## **SIK(SNF1LK)**

Product code: 02-131

Full-length human SIK [1-783(end) amino acids of accession number NP\_775490.2] was expressed as N-terminal GST-fusion protein (112 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: AMARA peptide
ATP ( $\mu$ M) Km app / Bin	: 47 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.8
IC50 at 1 mM ATP (nM)	: 1.0

---