

---

## QIK(SNF1LK2)

Product code 02-129

Full-length human QIK(SNF1LK2) [1-926(end) amino acids of accession number NP\_056006.1] was expressed as N-terminal GST-fusion protein (132 kDa) using baculovirus expression system. GST-QIK was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : AMARA peptide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 42 / 50

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 2.4

IC<sub>50</sub> at 1 mM ATP (nM) : 2.9

---