

---

---

## SIK3(QSK)

Product code: 02-130

Full-length human SIK3 [1-1369(end) amino acids of accession number NP\_001353615.1] was expressed as N-terminal GST-fusion protein (178 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: AMARA peptide
ATP ( $\mu$ M) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: 2.0

---

---