

---

## SGK3(SGKL)

Product code: 01-160

Truncated human SGK3 [119-496(end) amino acids and S486D of accession number NP\_037389.4] was co-expressed as N-terminal GST-fusion protein (68 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP\_002604.1] using baculovirus expression system. GST-SGK3 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: SGKtide
ATP ( $\mu$ M) Km app / Bin	: 17 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 42
IC50 at 1 mM ATP (nM)	: n.a.