

---

# SRPK1

Product code: 04-160

Full-length human SRPK1 [1-655(end) amino acids and V504 deletion of accession number NP\_003128.3] was expressed as N-terminal GST-fusion protein (101 kDa) using E. coli expression system.

Assay platform	: IMAP
Substrate	: SRPKtide
ATP ( $\mu$ M) Km app / Bin	: 200 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 85
IC50 at 1 mM ATP (nM)	: n.a.

---