

---

---

# BRSK2

Product code: 02-116

Full-length human BRSK2 [1-674(end) amino acids of accession number ABA17261.1] was expressed as N-terminal GST-fusion protein (102 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CHKtide
ATP ( $\mu$ M) Km app / Bin	: 31 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.31
IC50 at 1 mM ATP (nM)	: n.a.

