

---

---

# BRSK1

Product code: 02-115

Full-length human BRSK1 [1-778(end) amino acids of accession number NP\_115806.1] was expressed as N-terminal GST-fusion protein (112 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CHKtide
ATP ( $\mu$ M) Km app / Bin	: 30 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.27
IC50 at 1 mM ATP (nM)	: 0.57

