

---

# BRSK2

Product code: 02-116

Full-length human BRSK2 [1-674(end) amino acids of accession number NP\_001243558.1] was expressed as N-terminal GST-fusion protein (102 kDa) using baculovirus expression system.

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

---