
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. GST-BRSK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CHKtide

ATP (μ 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.
