
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 (μ 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 |

