

---

---

## BRK(PTK6)

Product code: 08-165

Full-length human BRK [2-451(end) amino acids of accession number NP\_005966.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP ( $\mu$ M) Km app / Bin	: 250 / 250
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 260
IC50 at 1 mM ATP (nM)	: 390

