

---

# BLK

Product code: 08-164

Full-length human BLK [1-505(end) amino acids of accession number NP\_001706.2] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 62 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.6
IC50 at 1 mM ATP (nM)	: 17

