

---

---

# PIP5K1B

Product code: 11-112

Full-length human PIP5K1B [1-540(end) amino acids of accession number NP\_003549.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system.

Assay platform : ADP-Glo

Substrate : PI(4)P

ATP ( $\mu$ M) Km app / Bin : 95 / 100

Metal : Mg

Reference compound : AG-183

IC50 at ATP Bin (nM) : 4300

IC50 at 1 mM ATP (nM) : n.a.