

---

---

# PIP5KL1

Product code: 11-114

Full-length human PIP5KL1 [1-394(end) amino acids of accession number NP\_001128691.1] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system.

Assay platform : ADP-Glo\_Lipid

Substrate : PI(4)P

ATP ( $\mu$ M) Km app / Bin : 1 / 1

Metal : Mg

Reference compound : AG-183

IC50 at ATP Bin (nM) : 2200

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

---

---