

---

---

# PIP4K2A

Product code: 11-115

Full-length human PIP4K2A [1-406(end) amino acids of accession number NP\_005019] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system.

Assay platform : ADP-Glo\_Lipid

Substrate : PI(5)P

ATP ( $\mu$ M) Km app / Bin : 20 / 25

Metal : Mg

Reference compound : AG-183

IC50 at ATP Bin (nM) : 7600

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

---

---