
PIP4K2B

Product code: 11-116

Full-length human PIP4K2B [1-416(end) amino acids of accession number NP_003550] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system.

Assay platform : ADP-Glo

Substrate : PI(5)P

ATP (μ M) Km app / Bin : 18 / 25

Metal : Mn

Reference compound : AG-183

IC50 at ATP Bin (nM) : 48000

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