
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. GST-PIP4K2B was purified by using glutathione sepharose chromatography.

Assay platform : ADP-Glo

Substrate : PI(5)P

ATP (μ M) K_{mapp} / B_{in} : 18 / 25

Metal : Mn

Reference compound : AG-183

IC₅₀ at ATP B_{in} (nM) : 48000

IC₅₀ at 1 mM ATP (nM) : n.a.
