
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

Substrate : PI(5)P

ATP (μ M) K_m app / B_{in} : 20 / 25

Metal : Mg

Reference compound : AG-183

IC₅₀ at ATP B_{in} (nM) : 7600

IC₅₀ at 1 mM ATP (nM) : n.a.