
PIP5K1A

Product code: 11-111

Full-length human PIP5K1A [1-549(end) amino acids of accession number NP_003548.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system.

Assay platform : ADP-Glo

Substrate : PI(4)P

ATP (μ M) K_m app / B_{in} : 28 / 25

Metal : Mg

Reference compound : AG-183

IC₅₀ at ATP B_{in} (nM) : 10000

IC₅₀ at 1 mM ATP (nM) : n.a.