
PIKFYVE(PIP5K3)

Product code: 11-118

Full-length human PIKFYVE [1-2098(end) amino acids and S696N, L932S, Q995L, T998S, S1033A and Q1183K of accession number NP_055855.2] was expressed as N-terminal GST-fusion protein (265 kDa) using baculovirus expression system.

Assay platform : ADP-Glo

Substrate : PI(3)P

ATP (μ M) Km app / Bin : 36 / 50

Metal : Mg

Reference compound : AG-183

IC50 at ATP Bin (nM) : 3900

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