
PDHK4(PDK4)

Product code: 10-125

Full-length human PDHK4 [1-411(end) amino acids of accession number NP_002603.1] was expressed as N-terminal GST-fusion protein (73 kDa) using E.coli expression system.

Assay platform : MSA
Substrate : PDHKtide

ATP (μ M) Km app / Bin : 19 / 25
Metal : Mg+K
Reference compound : VER-246608
IC50 at ATP Bin (nM) : 120
IC50 at 1 mM ATP (nM) : n.a.
