
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.

