
PDHK2(PDK2)

Product code: 10-140

Full-length human PDHK2 [1-407(end) amino acids of accession number NP_002602.2] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: PDHKtide
ATP (μ M) Km app / Bin	: 28 / 25
Metal	: Mg+K
Reference compound	: VER-246608
IC50 at ATP Bin (nM)	: 33
IC50 at 1 mM ATP (nM)	: n.a.

