
PHKG2

Product code: 02-153

Full-length human PHKG2 [1-406(end) amino acids of accession number NP_000285.1] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system.

| | |
|-------------------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : GS peptide |
| ATP (μ M) K_m app / B_{in} | : 8.1 / 10 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at ATP B_{in} (nM) | : 0.74 |
| IC50 at 1 mM ATP (nM) | : n.a. |

