
HCK

Product code: 08-169

Truncated human HCK [25-526(end) amino acids of accession number NP_002101.2] was expressed as N-terminal GST-fusion protein (84 kDa) using baculovirus expression system.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : Srctide |
| ATP (μ M) Km app / Bin | : 11 / 10 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 1.1 |
| IC50 at 1 mM ATP (nM) | : 22 |

