
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
