
CSK

Product code: 08-111

Full-length human CSK [1-450(end) amino acids of accession number NP_004374.1] was expressed as N-terminal GST-fusion protein (78 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Srctide

ATP (μ M) Km app / Bin : 4.8 / 5

Metal : Mg+Mn

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 53

IC50 at 1 mM ATP (nM) : 1500
