
PKAC β (PRKACB)

Product code: 01-128

Full-length human PKAC β [1-351(end) amino acids of accession number NP_002722.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : Kemptide

ATP (μ M) Km app / Bin : 4.7 / 5
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 1.0
IC50 at 1 mM ATP (nM) : 37
