
TAOK2

Product code: 07-133

Human TAOK2, catalytic domain [1-319 amino acid of accession number NP_004774.1] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : TAOKtide

ATP (μ M) Km app / Bin : 39 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 24
IC50 at 1 mM ATP (nM) : 100
