
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

