
MAPKAPK2

Product code: 02-142

Full-length human MAPKAPK2 [1-400(end) amino acids of accession number NP_116584.2] was co-expressed as N-terminal GST-fusion protein (73 kDa) with human His-tagged p38 β [1-364(end) amino acids of accession number NP_002742.3] and human His-tagged MAP2K6 [1-334(end) amino acids of accession number NP_002749.2] using baculovirus expression system.

Assay platform : MSA
Substrate : GS peptide

ATP (μ M) Km app / Bin : 3.6 / 5
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 80
IC50 at 1 mM ATP (nM) : 9300
