
MAPKAPK5

Product code: 02-144

Full-length human MAPKAPK5 [1-471(end) amino acids of accession number NP_003659.2] was co-expressed as N-terminal GST-fusion protein (81 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] using baculovirus expression system. GST-MAPKAPK5 was purified by using glutathione sepharose chromatography.

Assay platform : MSA
Substrate : GS peptide

ATP (μ M) Km app / Bin : 12 / 10
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 320
IC50 at 1 mM ATP (nM) : n.a.
