
MAPKAPK3

Product code: 02-143

Full-length human MAPKAPK3 [1-382(end) amino acids of accession number NP_004626.1] was co-expressed as N-terminal GST-fusion protein (70 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. GST-MAPKAPK3 was purified by using glutathione sepharose chromatography.

Assay platform : MSA
Substrate : GS peptide

ATP (μ M) Km app / Bin : 13 / 10
Metal : Mg
Reference compound : K252a
IC50 at ATP Bin (nM) : 410
IC50 at 1 mM ATP (nM) : n.a.
