
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 and Ni-NTA chromatography.

Assay platform : Mobility Shift Assay

Substrate : GS peptide

ATP (μ M) K_{mapp} / B_{in} : 13 / 10

Metal : Mg

Reference compound : K252a

IC₅₀ at ATP B_{in} (nM) : 410

IC₅₀ at 1 mM ATP (nM) : n.a.