
MAP3K2

Product code: 07-104

Human MAP3K2, catalytic domain [336-619(end) amino acids of accession number NP_006600.3] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : MAP2K4/MAP2K7
Cascade Assay *
ATP (μ M) : 1000
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 45

* (MAP2K4/MAP2K7)/JNK2/Modified Erktide
