
MINK(MINK1)

Product code: 07-139

Human MINK, catalytic domain [1-314 amino acids of accession number NP_056531.1] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : Modified Erktide

ATP (μ M) Km app / Bin : 16 / 50

Metal : Mg

Reference compound : K252a

IC50 at ATP Bin (nM) : 5.5

IC50 at 1 mM ATP (nM) : 4.7
