
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
