
p38 β (MAPK11)

Product code: 04-153

Full-length human p38 β [1-364(end) amino acids of accession number NP_002742.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system.

Assay platform	: MSA
Substrate	: Modified Erktide
ATP (μ M) Km app / Bin	: 63 / 75
Metal	: Mg
Reference compound	: SB202190
IC50 at ATP Bin (nM)	: 16
IC50 at 1 mM ATP (nM)	: 110
