
MARK2

Product code 02-121

Full-length human MARK2 [1-745(end) amino acids of accession number NP_059672.2] was expressed as N-terminal GST-fusion protein (110 kDa) using baculovirus expression system. GST-MARK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CHKtide

ATP (μ M) K_{mapp} / B_{in} : 8.8 / 10

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 0.12

IC₅₀ at 1 mM ATP (nM) : n.a.
