
MARK1

Product code 02-120

Full-length human MARK1 [1-795(end) amino acids of accession number AAF72103.1] was expressed as N-terminal GST-fusion protein (116 kDa) using baculovirus expression system. GST-MARK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CHKtide

ATP (μ M) K_{mapp} / B_{in} : 8 / 10

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 0.17

IC₅₀ at 1 mM ATP (nM) : n.a.
