
ITK

Product code 08-181

Full-length human ITK [2-620(end) amino acids of accession number NP_005537.3] was expressed as N-terminal GST-fusion protein (99 kDa) using baculovirus expression system. GST-ITK was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP (μ M) K_m app / B_{in} : 6.1 / 10

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 3.2

IC₅₀ at 1 mM ATP (nM) : 200
