
SRC

Product code 08-173

Full-length human SRC [1-536(end) amino acids of accession number NP_005408.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-SRC was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Srctide

ATP (μ M) K_{mapp} / B_{in} : 31 / 50

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 5.3

IC₅₀ at 1 mM ATP (nM) : 33
