
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.

Assay platform : MSA

Substrate : Srctide

ATP (μ M) Km app / Bin : 31 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 5.3

IC50 at 1 mM ATP (nM) : 33
