
ROCK1

Product code: 01-109

Human ROCK1, catalytic domain [1-477 amino acids of accession number NP_005397.1] was expressed as N-terminal GST-fusion protein (82 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : LIMKtide

ATP (μ M) K_m app / B_{in} : 3.1 / 5
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP B_{in} (nM) : 1.6
IC50 at 1 mM ATP (nM) : 73
