
ROCK2

Product code: 01-110

Human ROCK2, catalytic domain [1-553 amino acids of accession number NP_004841.2] was expressed as N-terminal GST-fusion protein (91 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: LIMKtide
ATP (μ M) K_m app / B_{in}	: 7.4 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP B_{in} (nM)	: 0.92
IC50 at 1 mM ATP (nM)	: 21

