
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
