
RAF1

Product code: 09-125

Human RAF1, catalytic domain [306-648(end) amino acids and Y340D and Y341D of accession number NP_002871.1] was expressed as N-terminal GST-fusion protein (66 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : MAP2K1
Cascade Assay *
ATP (μ M) : 1000
Metal : Mg
Reference compound : ZM336372
IC50 at 1 mM ATP (nM) : >10000

* MAP2K1/Erk2/Modified Erktide
