
BRAF

Product code: 09-122

Human BRAF, catalytic domain [433-726 amino acid of accession number NP_004324.2] was expressed as N-terminal GST-fusion protein (60 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
