
BRAF[V600E]

Product code: 09-144

Human BRAF, catalytic domain [433-726 amino acids and V600E 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)	:	662

* MAP2K1/Erk2/Modified Erktide
