

---

# 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 (μM)	: 1000
Metal	: Mg
Reference compound	: ZM336372
IC50 at 1 mM ATP (nM)	: >10000

\* MAP2K1/Erk2/Modified Erktide

---