





## RafRBD (c-Raf)

Ras Binding Domain of Raf, c-Raf 1 Kinase human, recombinant, *E. coli* 

Cat. No.	Amount
PR-305	50 µg

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 9.4 kDa (50 - 132 amino acids)

Accession number: X03484 / P04049

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 32 mM Tris-HCl pH 7.6, 200 mM NaCl, 5 mM  $\rm MgCl_2$  and 5 mM DTT)

## Description:

Raf is a member of the serine/threonine proteine kinase family involved in regulation of cell growth and differentiation and is the most important effector of Ras. Full-length Raf is composed of three conserved regions responsible for interaction with Ras, for phosphorylation and for catalytic activity. RafRBD (Ras binding domain, amino acids 50-132) mediates interaction with membrane-anchored Ras necessary for activation of the kinase activity of Raf.

## Selected References:

Eing *et al.* (2002) Quantification of the Raf-C1 interaction with solidsupported bilayers. *ChemBioChem* **3**:190.

Emerson *et al.* (1996) Structure of Ras-binding domain of c-Raf-1 as determined by NMR spectroscopy and identification of the region that interacts with Ras. *Drug Des. Discov.* **13**:83.

Nasser *et al.* (1995) The 2.2 crystal structure of the Ras-binding domain of the serine/threonine kinase c-Raf1 in complex with Rap1A and a GTP analogue. *Nature* **375**:554.

Bahk *et al.* (2008) A Cross-talk between oncogenic Ras and tumor suppressor PTEN through FAK Tyr861 phosphorylation in NIH/3T3 mouse embryonic fibroblasts. *Biochemical and Biophysical Research Communications* **377** (4):1199-1204.

