

## RafRBD (c-Raf)

human, recombinant, *E. coli*

Cat. No.	Amount
PR-305	50 µg

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -80°C

### Avoid freeze / thaw cycles

### Form

Liquid. Supplied in 32 mM Tris-HCl pH 7.6, 200 mM NaCl, 5 mM MgCl<sub>2</sub> and 5 mM DTT.

### Molecular Weight

9.4 kDa (amino acids 50 - 132)

### Purity

> 95% by SDS-PAGE

### Description

Raf is a member of the serine/threonine protein 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.

### Associated products available from Jena Bioscience

**Fluorescent ATPs** (such as MANT, ANT, TNP, or ATP labeled with a dye of your choice)

**Non-hydrolyzable ATP-analogs** (such as ATP-γS, AppCp, AppNHp, NPE-caged-ATP...)

**Recombinant G-protein signaling reagents** (including GAPs, GEFs, small GTPases, heterotrimeric G-proteins, and many more...)

**Kinase Substrate Peptides**

For further information please visit [www.jenabioscience.com](http://www.jenabioscience.com).