





RafRBD^{GST} (C-Raf)

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

Cat. No.	Amount
PR-366	300 µg

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: thaw quickly/avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 42 kDa

Accession number: X03484 / P04049

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in PBS, pH 7.4)

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. The GST-Tag facilitates the protein's application in typical GST pull-down assays. The protein is GST-tagged on its N-terminus. There is a 7 AA linker between GST and RBD.

Reommended thawing and freezing :

Thaw the protein quickly at approx. 30 °C in a water bath and refreeze aliquots fast on dry ice.

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.

