

RalGDS-RBD^{GST}

Ras Binding Domain of the Ral Guanine Nucleotide Dissociation Stimulator
human, recombinant, *E. coli*

Cat. No.	Amount
PR-365	0.25 mg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 30 mM Tris-HCl pH 7.5, 100 mM NaCl, 5 mM MgCl₂ and 5 mM DTE.

Purity

≥90% by SDS-PAGE

Description

Ras-binding domain (RBD) of the human RalGDS, comprises amino acids 1 - 97 of wild type.

RalGDS (Ral-guanine nucleotide dissociation stimulator) is the guanine nucleotide exchange factor (GEF) for the Ras-related protein Ral. The GEF-activity of RalGDS is stimulated by binding of Ras.

The GST-tag facilitates the protein's application in typical GST pull down assays.

Selected References:

Vetter *et al.* (1999) Structural and biochemical analysis of Ras effector signaling via RalGDS. *FEBS Letters* **451**:175.

Murphy *et al.* (2002) Involvement of phosphatidylinositol 3-kinase, but not RalGDS, in TC21/R-Ras2-mediated transformation. *J. Biol. Chem.* **277**:9966.