

## RalGDS-RBD<sup>GST</sup>

(Ras Binding Domain of the Ral Guanine Nucleotide Dissociation Stimulator)

Human, Recombinant, *E. coli*

Cat. No.	Amount
PR-365	0.25 mg

Liquid. Supplied as a 1 mg/ml solution in 30 mM Tris-HCl, pH 7.5, 100 mM NaCl, 5 mM MgCl<sub>2</sub>, and 5 mM DTE.

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.

AVOID FREEZE/THAW CYCLES!

**For in vitro use only!**

**Purity:** > 90% by SDS-PAGE

**Store:** -80 °C

### 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.