

**RhotekinRBD^{GST}**

Rhotekin Ras Binding Domain, from the Japanese human, recombinant, *E. coli*

Cat. No.	Amount
PR-367	50 µg

For in vitro use only!

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Accession number: NM_033046

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 50 mM Tris-HCl pH 7.5, 100 mM NaCl and 2 mM DTE)

Description:

Rhotekin (from the Japanese teki, meaning target) is an effector of the Rho-GTPases that was previously identified in mice. GST-RhotekinRBD is a glutathione S-transferase (GST)-fusion protein of the Rho-binding domain (RBD) of Rhotekin. The N-terminal peptide of Rhotekin contains the amino acids 1-89 comprising the Rho-binding domain between amino acids 7 and 89. The GST-Tag facilitates the protein's application in typical GST pull down assays.

Selected References:

Reynaud *et al.* (2000) The PDZ protein TIP-1 interacts with the Rho effector Rhotekin and is involved in Rho signaling to the serum response element. *J. Biol. Chem.* **275**:33962.

Reid *et al.* (1996) Rhotekin, a new putative target for Rho bearing homology to the serine/threonine kinase, PKN, and rhodophilin in the Rho-binding domain. *J. Biol. Chem.* **271**:13556.