

## RhotekinRBD<sup>GST</sup>

(Rhotekin Ras Binding Domain, from the Japanese Human, Recombinant, *E. coli*)

Cat. No.	Amount
PR-367	50 µg

Liquid. Supplied as a 1 mg/ml solution in 50 mM Tris-HCl, pH 7.5, 100 mM NaCl, and 2 mM DTE.

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.

AVOID FREEZE/THAW CYCLES!

**For in vitro use only!**

**Purity:** > 95% by SDS-PAGE

**Store:** -80 °C

### 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.