

## AF6-RBD<sup>GST</sup> (residues 1-141)

(GTP-H-Ras binding cytoplasmic protein, Afadin - Ras-binding domain)

Human, Recombinant, *E. coli*

Cat. No.	Amount
PR-398	50 µg

Liquid. Supplied in 25 mM Hepes, pH 7.5, 50 mM NaCl, 1 mM EDTA, and 50% glycerol.

The amino terminus of Af6 was identified as the Ras-binding site. Additionally to the Ras-binding Af6 also binds Rap. ZO-1, a protein involved in the formation of tight junctions, also binds to Af6 close to the amino terminus thereby competing with Ras binding. These data suggest a participation of Af6 in the regulation of cell-cell contacts via a Ras-modulated interaction with ZO-1. In contrast to the well characterized effectors Raf, RalGEF, and PI(3)kinase, Af6 is a protein with no enzymatic function. It seems to serve as a scaffolding component for protein complexes. This would represent a new type of Ras effector because Ras binding in this case does not lead to the activation of the enzymatic activity of the effector.

AVOID FREEZE/THAW CYCLES!

**For in vitro use only!**

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

**Store:** -20 °C

### Selected References:

Linnemann *et al.* (1999) Thermodynamic and Kinetic Characterization of the Interaction between the Ras Binding Domain of AF6 and Members of the Ras Subfamily. *J. Biol. Chem.* **274**:13556.

Steiner *et al.* (2000) Sequence-specific resonance assignment of the Ras-binding domain of AF6. *J. Biomol. NMR.* **18**:73.

### Associated products available from Jena Bioscience

**Non-hydrolyzable GTP-analogs** (such as GTPγS, GppCp, GppNHp, NPE-caged-GTP...)

**Fluorescent GTPs** (such as MANT, ANT, TNP, or GTP labeled with a dye of your choice)

**Recombinant G-protein signaling reagents** including GAPs, GEFs, heterotrimeric G-proteins, and many more...

**For detailed information please view the sections on**

<http://www.jenabioscience.com>