

AF6-RBD^{GST} (residues 1-141)

GTP-H-Ras binding cytoplasmic protein, Afadin - Ras-binding domain
human, recombinant, *E. coli*

Cat. No.	Amount
PR-398	50 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

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

Purity

> 95% by SDS-PAGE

Description

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.

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

Fluorescent GTPs (such as MANT, ANT, TNP, or GTP

labeled with a dye of your choice)

Non-hydrolyzable GTP-analogs (such as GTP γ S, GppCp, GppNHp, NPE-caged-GTP...)

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

For further information please visit www.jenabioscience.com.