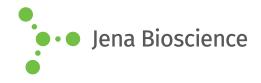
## **DATA SHEET**





## 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

For general laboratory use.

**Shipping:** shipped on gel packs **Storage Conditions:** store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

**Accession number:** NM\_005936

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 25 mM HEPES pH 7.5, 50 mM NaCl, 1 mM

EDTA and 50% glycerol)

## **Description:**

The amino terminus of Af6 was identified as the Rasbinding 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.