

## Recombinant ADAR1 (1-295) protein

**Catalog No:** 82140, 82740

**Expressed In:** *E. coli*

**Quantity:** 100, 1000 µg

**Source:** Human

**Buffer Contents:** Recombinant ADAR1 (1-295) protein is supplied in 25 mM Tris pH 8.0, 300 mM NaCl, 20% glycerol, 0.5 mM TCEP.

**Background:** ADAR1 (Double-stranded RNA-specific adenosine deaminase), Catalyzes the hydrolytic deamination of adenosine to inosine in double-stranded RNA (dsRNA) referred to as A-to-I RNA editing. This may affect gene expression and function in a number of ways that include mRNA translation by changing codons and hence the amino acid sequence of proteins since the translational machinery read the inosine as a guanosine; pre-mRNA splicing by altering splice site recognition sequences; RNA stability by changing sequences involved in nuclease recognition; genetic stability in the case of RNA virus genomes by changing sequences during viral RNA replication; and RNA structure-dependent activities such as microRNA production or targeting or protein-RNA interactions. Can edit both viral and cellular RNAs and can edit RNAs at multiple sites (hyper-editing) or at specific sites (site-specific editing).

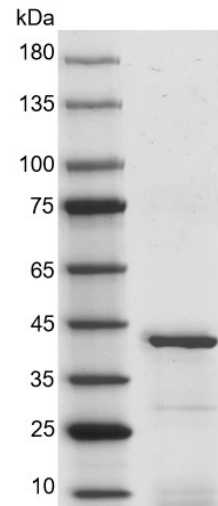
**Protein Details:** Recombinant ADAR1 (1-295) protein that includes amino acids 1-295 of human ADAR1 protein (accession number NP\_001102.3) was expressed in *E. coli* cells with an N-terminal 6xHis and Flag tag. The molecular weight of the protein is 35.88 kDa.

**Application Notes:** This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for this product is shown.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

### ADAR1 (1-295)



### Recombinant ADAR1 (1-295) protein

4-20% SDS-PAGE Coomassie staining

MW: 35.88 kDa,

Purity: ≥ 95%