

# PRODUCT INFORMATION



## ADAR2 Long Isoform (human, recombinant)

Item No. 42937

### Overview and Properties

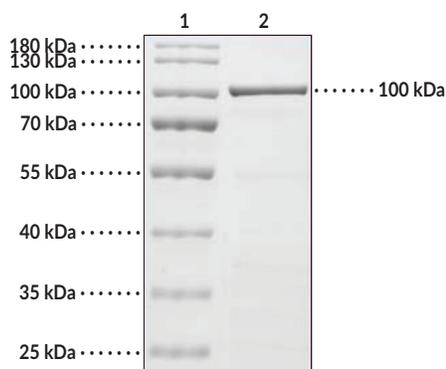
**Synonyms:** ADARB1, Adenosine Deaminase RNA-specific B1, DRADA2B  
**Source:** Recombinant human N-terminal DYKDDDDK-tagged ADAR2L expressed in insect cells  
**Amino Acids:** 2-741 (full length)  
**Uniprot No.:** P78563  
**Molecular Weight:** ~85 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** ≥85% estimated by SDS-PAGE  
**Supplied in:** 50 mM Tris, pH 7.5, with 150 mM sodium chloride, 1 mM EDTA, 0.5 mg/ml DYKDDDDK-tag, 1 mM DTT, and 10% glycerol

### Protein

**Concentration:** *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: ADAR2 Long Isoform

**SDS-PAGE Analysis of ADAR2 Long Isoform.** This protein has a calculated molecular weight of approximately 85 kDa. It has an apparent molecular weight of approximately 100 kDa by SDS-PAGE.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/12/2025

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Adenosine deaminase acting on RNA 2 (ADAR2) is an aminohydrolase and member of the ADAR protein family.<sup>1,2</sup> It is composed of an N-terminal nuclear export (NES) region, two RNA-binding domains, and a C-terminal deaminase domain.<sup>2</sup> ADAR2 is widely expressed and its two predominant isoforms, ADAR2 short (ADAR2S) and ADAR2 long (ADAR2L), form homodimers and localize primarily to the nucleus.<sup>1,3</sup> ADAR2L performs site-specific editing of dsRNA by converting adenosine to inosine, a post-transcriptional modification known as A-to-I editing that predominantly targets RNA encoding neuroreceptors and ion channels. ADAR2 has a variety of RNA substrates, including *Gria2*, which encodes the AMPA receptor GluA2 subunit, also known as GluR-B, and the adenosine at the Q/R site is edited by ADAR2 in nearly all transcripts, resulting in a codon change from CAG (glutamine) to CIG (arginine).<sup>3-5</sup> This edit reduces calcium permeability of the AMPA receptor and reduces trafficking of the receptor to the postsynaptic membrane.<sup>3,5</sup> Mice lacking *AdarB1*, the gene encoding *Adar2*, have reduced *Gria2* pre-mRNA editing, develop early-onset epilepsy, and are viable only up to three weeks.<sup>5</sup> Biallelic *ADARB1* variants are associated with intractable epilepsy, microcephaly, and intellectual disability.<sup>3</sup> *ADARB1* expression is decreased in ovarian cancer tumors and correlates with poor prognosis and shorter progression-free survival.<sup>6</sup> Cayman's ADAR2L (human, recombinant) protein has a calculated molecular weight of approximately 85 kDa. By SDS-PAGE, the apparent molecular mass of the protein is approximately 100 kDa.

## References

---

1. Zinshteyn, B. and Nishikura, K. Adenosine-to-inosine RNA editing. *Wiley Interdiscip. Rev. Syst. Biol. Med.* **1(2)**, 202-209 (2009).
2. Rehwinkel, J. and Mehdipour, P. ADAR1: From basic mechanisms to inhibitors. *Trends Cell Biol.* **35(1)**, 59-73 (2025).
3. Tan, T.Y., Sedmík, J., Fitzgerald, M.P., et al. Bi-allelic *ADARB1* variants associated with microcephaly, intellectual disability, and seizures. *Am. J. Hum. Genet.* **106(4)**, 467-483 (2020).
4. Macbeth, M.R., Schubert, H.L., Vandemark, A.P., et al. Inositol hexakisphosphate is bound in the ADAR2 core and required for RNA editing. *Science* **309(5740)**, 1534-1539 (2005).
5. Higuchi, M., Maas, S., Single, F.N., et al. Point mutation in an AMPA receptor gene rescues lethality in mice deficient in the RNA-editing enzyme ADAR2. *Nature* **406(6791)**, 78-81 (2000).
6. Zhu, W., Xu, Z., Huang, M., et al. Downregulated *ADARB1* facilitates cell proliferation, invasion and has effect on the immune regulation in ovarian cancer. *Front. Bioeng. Biotechnol.* **9:792911**, (2021).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM