

# PRODUCT INFORMATION



## YTHDF2 (human, recombinant)

Item No. 26345

### Overview and Properties

**Synonyms:** CAHL, CLL-associated Antigen KW-14, HGRG8, High-Glucose-Regulated Protein 8, Renal Carcinoma Antigen NY-REN-2, YTH Domain-containing Family Member 2, YTH N<sup>6</sup>-Methyladenosine RNA Binding Protein 2 (358-502; YTH domain)

**Source:** Recombinant N-terminal GST-tagged YTHDF2 purified from *E. coli*

**Amino Acids:** 358-502

**Uniprot No.:** Q9Y5A9

**Molecular Weight:** 43.24 kDa

**Storage:** -80°C (as supplied)

**Stability:** ≥2 years

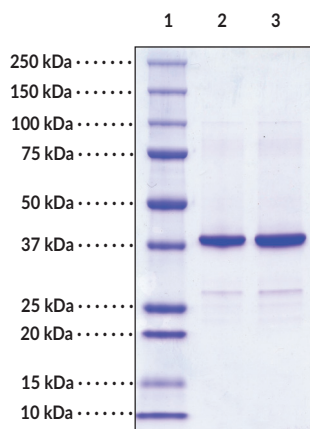
**Purity:** *batch specific* (≥90% estimated by SDS-PAGE)

**Supplied in:** 50 mM HEPES, pH 7.5, with 150 mM sodium chloride 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: YTHDF2 (2 µg)  
Lane 3: YTHDF2 (4 µg)

Representative gel image shown; actual purity may vary between each batch.

#### 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, 12/03/2025

#### 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

---

YTH-Domain family member 2 (YTHDF2) is a reader protein that binds to N<sup>6</sup>-methyladenosine (m<sup>6</sup>A) to regulate mRNA degradation.<sup>1,2</sup> YTHDF2 binds to m<sup>6</sup>A sites in the 3'-UTR of mRNA through its C-terminal YTH domain and recruits the CCR4-NOT deadenylase complex via its N-terminus to accelerate mRNA localization to decay sites, such as P-bodies, and promote subsequent degradation.<sup>1-3</sup> *Ythdf2* knockout results in a build-up of maternal mRNA in oocytes that induces female-specific infertility in mice.<sup>4</sup> Conditional knockout of *Ythdf2* increases the number of hematopoietic stem cells (HSCs) without shifting lineage differentiation or inducing malignancies.<sup>5</sup> It also increases *ex vivo* expansion of human umbilical cord HSCs. This product is the C-terminus of YTHDF2 corresponding to residues 358-502 of the human full-length sequence.

## References

---

1. Du, H., Zhao, Y., He, J., *et al.* YTHDF2 destabilizes m<sup>6</sup>A-containing RNA through direct recruitment of the CCR4-NOT deadenylase complex. *Nat. Commun.* **7:12626** (2016).
2. Zhong, L., Liao, D., Zhang, M., *et al.* YTHDF2 suppresses cell proliferation and growth via destabilizing the EGFR mRNA in hepatocellular carcinoma. *Cancer Lett.* **442**, 252-261 (2019).
3. Wang, X., Lu, Z., Gomez, A., *et al.* N<sup>6</sup>-Methyladenosine-dependent regulation of messenger RNA stability. *Nature* **505(7481)**, 117-120 (2014).
4. Ivanova, I., Much, C., Di Giacomo, M., *et al.* The RNA m<sup>6</sup>A reader YTHDF2 is essential for the post-transcriptional regulation of the maternal transcriptome and oocyte competence. *Mol. Cell* **67(6)**, 1059-1067 (2017).
5. Li, Z., Qian, P., Shao, W., *et al.* Suppression of m<sup>6</sup>A reader *Ythdf2* promotes hematopoietic stem cell expansion. *Cell Res.* **28(9)**, 904-917 (2018).