

## Recombinant YTHDC2 protein

---

**Catalog No:** 81198

**Lot No:** 25018001

**Expressed In:** Baculovirus

**Quantity:** 20 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant YTHDC2 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

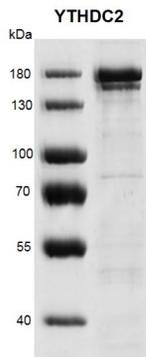
**Background:** YTHDC2 (YTH Domain Containing 2), is a member of the DEAH (Asp-Glu-Ala-His) subfamily of proteins, part of the DEAD (Asp-Glu-Ala-Asp) box family of RNA helicases. It specifically recognizes and binds to N6-methyladenosine (m6A), a common modified RNA nucleotide that is enriched in the stop codons and 3' UTRs of eukaryotic mRNAs and some non-coding RNAs and plays a role in the efficiency of mRNA splicing, processing and stability. YTHDC2

may be associated with susceptibility to pancreatic cancer in human patients, and knockdown of this gene resulted in reduced proliferation in a human liver cancer cell line.

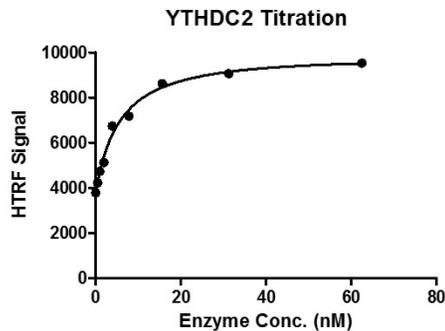
**Protein Details:** Recombinant YTHDC2 protein was expressed in a baculovirus expression system as the full length protein (accession number NP\_073739.3) with an N-terminal Flag tag. The molecular weight of the protein is 161.5 kDa.

**Application Notes:** Recombinant YTHDC2 protein is suitable for use in the study of binding assay, inhibitor screening, and selectivity profiling.

**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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

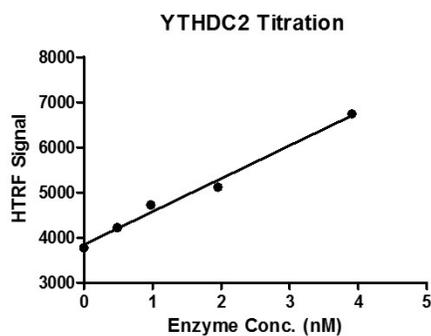


**Recombinant YTHDC2 protein gel**  
 8% SDS-PAGE Coomassie staining  
 MW: 161.5 kDa  
 Purity: >87%



### Recombinant YTHDC2 binding assay

3  $\mu$ M biotinylated ssDNA m6A oligos (GTTGG/m6A/CTT) were incubated with different concentrations of YTHDC2 protein in a 10  $\mu$ l system containing 50 mM HEPES-NaOH pH 7.4, and 0.1% BSA for 1 hour, then 10  $\mu$ l FLAG antibody and SA-XL665 mixture (each 1:100 dilution reaction buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.



### Recombinant YTHDC2 binding assay

3  $\mu$ M biotinylated ssDNA m6A oligos (GTTGG/m6A/CTT) were incubated with different concentrations of YTHDC2 protein in a 10  $\mu$ l system containing 50 mM HEPES-NaOH pH 7.4, and 0.1% BSA for 1 hour, then 10  $\mu$ l FLAG antibody and SA-XL665 mixture (each 1:100 dilution reaction buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.