

YTHDF2 Polyclonal Antibody

(Catalog # A68190)

Background

Specifically recognizes and binds N6-methyladenosine (m6A)-containing RNAs, and regulates mRNA stability (PubMed:24284625, PubMed:26046440). M6A is a modification present at internal sites of mRNAs and some non-coding RNAs and plays a role in the efficiency of mRNA splicing, processing and stability (PubMed:22575960, PubMed:24284625, PubMed:25412658, PubMed:25412661). Acts as a regulator of mRNA stability: binding to m6A-containing mRNAs results in the localization to mRNA decay sites, such as processing bodies (P-bodies), leading to mRNA degradation (PubMed:24284625, PubMed:26046440).

Description

YTHDF2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q9Y5A9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human YTH domain-containing family protein 2 protein (201-500AA)

Storage

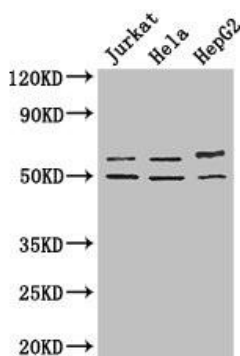
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

YTH domain-containing family protein 2, CLL-associated antigen KW-141, YTHDF2

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:2000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

Positive WB detected in: Jurkat whole cell lysate, HeLa whole cell lysate, HepG2 whole cell lysate

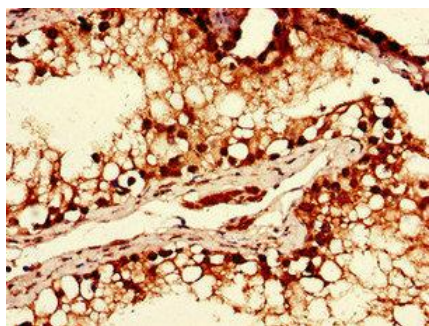
All lanes: YTHDF2 antibody at 2.7ug/ml

Secondary

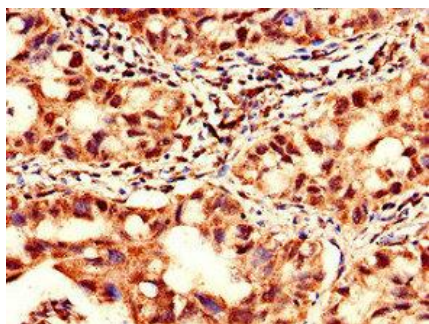
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 63, 57 kDa

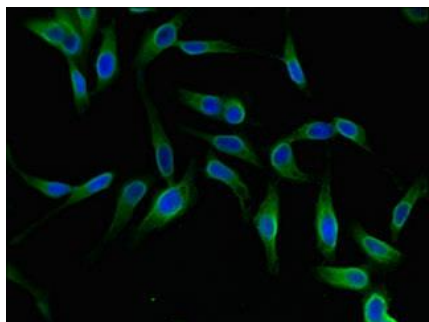
Observed band size: 63, 50 kDa



Immunohistochemistry of paraffin-embedded human testis tissue using YTHDF2 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human lung cancer using YTHDF2 Antibody at dilution of 1:100



Immunofluorescent analysis of HeLa cells using YTHDF2 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)