
HNRNPA1 Polyclonal Antibody

(Catalog # A73393)

Background

This gene encodes a member of a family of ubiquitously expressed heterogeneous nuclear ribonucleoproteins (hnRNPs), which are RNA-binding proteins that associate with pre-mRNAs in the nucleus and influence pre-mRNA processing, as well as other aspects of mRNA metabolism and transport. The protein encoded by this gene is one of the most abundant core proteins of hnRNP complexes and plays a key role in the regulation of alternative splicing. Mutations in this gene have been observed in individuals with amyotrophic lateral sclerosis 20. Multiple alternatively spliced transcript variants have been found. There are numerous pseudogenes of this gene distributed throughout the genome.

Description

HNRNPA1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P09651

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-372 of human HNRNPA1 (NP_112420.1).

Storage

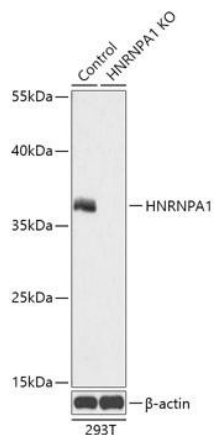
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

HNRNPA1; ALS19; ALS20; HNRPA1; HNRPA1L3; IBMPFD3; UP 1; hnRNP A1; hnRNP-A1

Application

WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200



Western blot analysis of extracts from normal (control) and HNRNPA1 knockout (KO) 293T cells, using HNRNPA1 Polyclonal Antibody at 1:200 dilution.

Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 10s.