

Inosine antibody

Cat. No. C15200251

Type: Monoclonal

Isotype: IgG1

Source: Mouse

Lot: 002

Size: 50 µg

Concentration: 2 µg/µl

Specificity: Human, other (wide range): positive.

Purity: Protein A purified monoclonal antibody.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Storage buffer: PBS containing 0.05% sodium azide.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Monoclonal antibody raised in mouse against Inosine (I) conjugated to BSA.

Applications

Applications	Suggested dilution	References
RIP*	1 µg per IP	Fig 1

*Please note that the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-10 µg per IP.

Target description

The formation of Inosine by deamination of adenosine is probably one of the most common post-translational modifications in RNA. It is an essential modification introduced by specialized enzymes in a highly regulated manner to generate transcriptome diversity and deficiencies of the enzymes involved in this conversion lead to a variety in diseases including cancer, viral infections and neurological and psychiatric disorders. Only a small part of the Inosine conversions occurs in coding sequences; the vast majority is present in noncoding sequences such as microRNAs, tRNAs, and introns and 3' untranslated regions of messenger RNAs, which play important roles in the RNA-mediated regulation of gene expression.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue du Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode LLC. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: January, 2023

Results

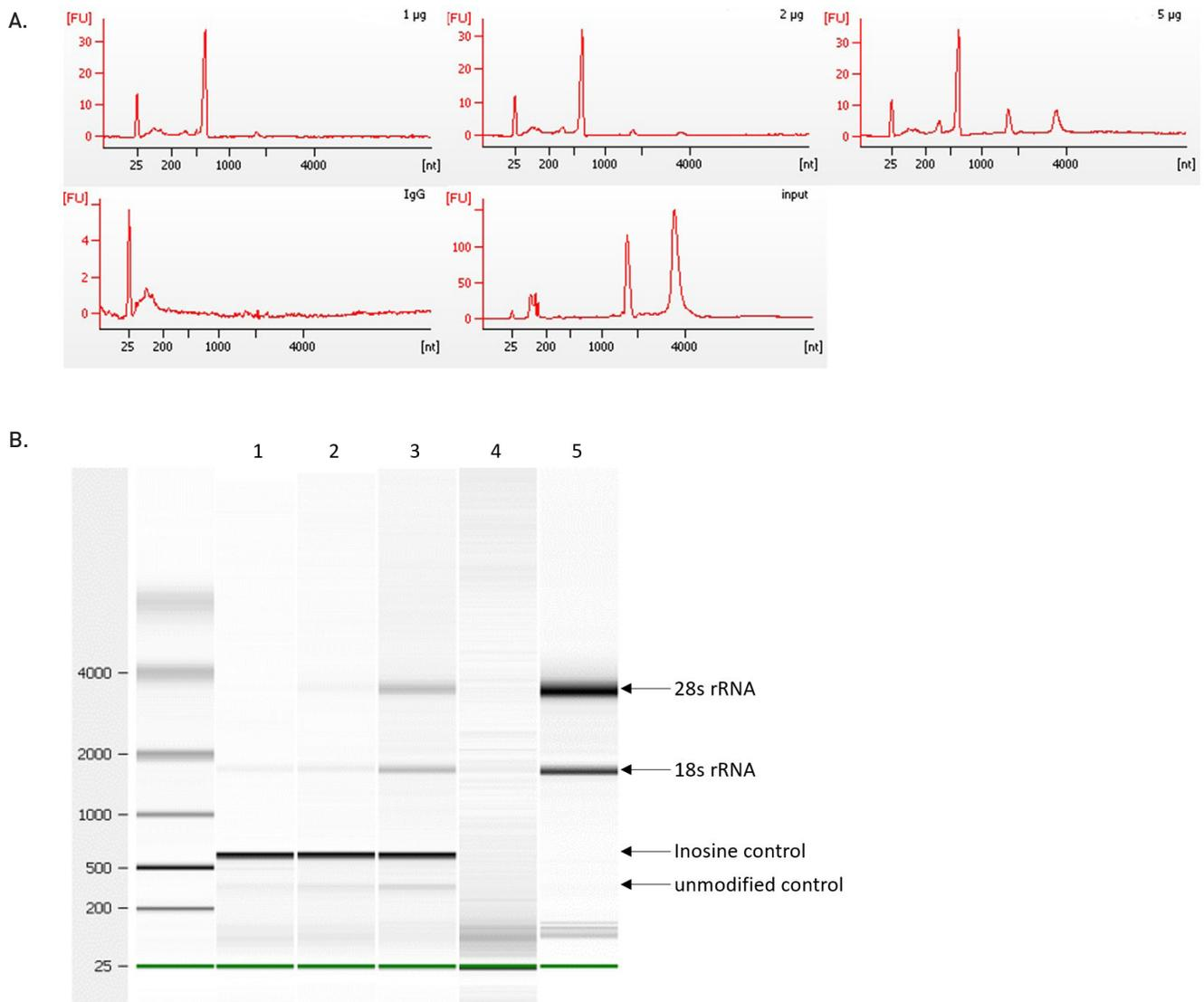


Figure 1. RNA immunoprecipitation using the Diagenode monoclonal antibody directed against Inosine

IP was performed with the Diagenode antibody against Inosine (cat. No. C15200251) on 40 µg total RNA from human HeLa cells, spiked with an in vitro produced RNA molecule containing Inosine nucleotides as well as an unmodified control RNA (40 ng each). A titration of the antibody of 1, 2 and 5 µg was analyzed. IgG (2 µg/IP) was used as negative control.

Figure 1A The immunoprecipitated RNA was subsequently analysed on a Bioanalyzer. The peak at ~500 bp corresponds to the Inosine spike in, whereas the negative control spike in (expected size ~300 bp) is not captured by the antibody.

Figure 1B shows the gel image for the Inosine antibody (lane 1, 2 and 3), the IgG negative control (lane 4) and the input (lane 5). The marker (in bp) is shown on the left, the position of the 28s and 18s ribosomal RNA and of both spike in controls is indicated on the right.