

DROSHA polyclonal antibody

Other names: RN3, RNASEN, RNASE3L, RNase III, p241, RANSE3L, ETOHI2

Cat. No. C15310246 (CS-PA026-100)

Type: Polyclonal

Source: Rabbit

Lot #: A1171-004

Size: 100 µl

Concentration: not determined

Specificity: Mouse: positive / Other species: not tested

Purity: Whole antiserum from rabbit containing 0.05% azide.

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

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

Description: Polyclonal antibody raised in rabbit against mouse DROSHA (drosha, ribonuclease type III) using two KLH-conjugated synthetic peptides containing a sequence from the central and the C-terminal region of the protein, respectively.

Applications

	Suggested dilution	Results
ELISA	1:500	Fig 1

*The optimal dilution for other applications should be determined by the end user. For WB we suggest starting with a 1:1,000 dilution

Target description

DROSHA (UniProtKB/Swiss-Prot entry Q9NRR4) (RNase III) is a double-stranded RNA-specific endoribonuclease that is involved in the initial step of microRNA (miRNA) biogenesis. DROSHA is a component of the microprocessor complex that is required to process primary miRNA transcripts (pri-miRNAs) to release precursor miRNA (pre-miRNA) in the nucleus. Within this complex, DROSHA cleaves the 3' and 5' strands of a stem-loop in pri-miRNAs (processing center 11 bp from the dsRNA-ssRNA junction) to release hairpin-shaped pre-miRNAs that are subsequently cut by the cytoplasmic DICER to generate mature miRNAs. Involved also in pre-rRNA processing. Cleaves double-strand RNA and does not cleave single-strand RNA. Involved in the formation of GW bodies

Results

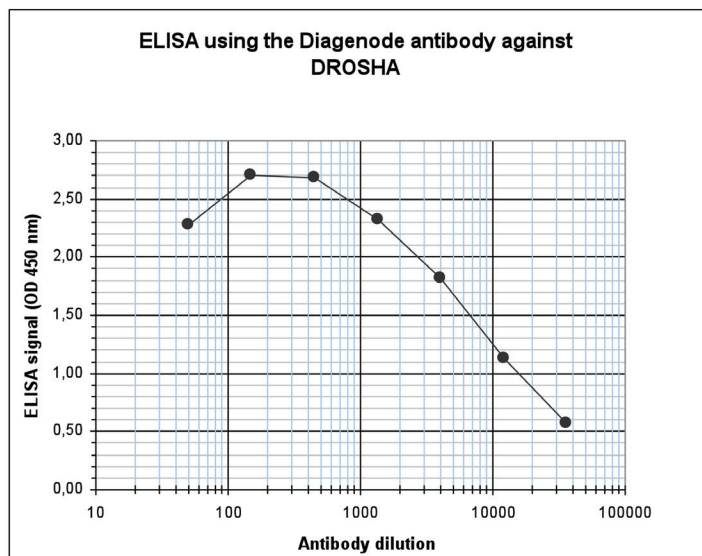


Figure 1. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against mouse DROSHA (cat. No. CS-PA026-100). The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:8,100.

Diagenode sa. BELGIUM | EUROPE

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

Diagenode Inc. 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: September 5, 2012