

hnRNP A/B polyclonal antibody - Classic

Cat. No. C15410263

Type: Polyclonal

Source: Rabbit

Lot #: 39645

Size: 25 µl/100 µl

Concentration: 1 µg/µl

Specificity: Human.

Purity: Purified by peptide affinity chromatography.

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.

Applications

	Suggested dilution*	Results
ICC/IF	1:100-1:1,000	Figure 3
IHC	1:100-1:1,000	Figure 1
Western blot	1:500-1:3,000	Figure 2

*Optimal dilutions/concentrations should be determined by the researcher.

Results

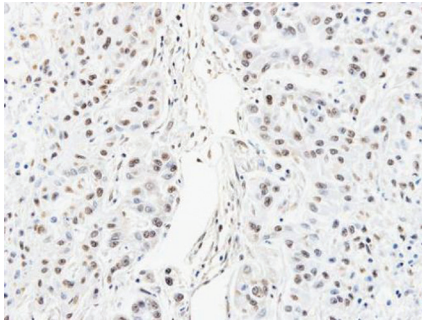


Figure 1. IHC

Immunohistochemical analysis of paraffin-embedded A549 xenograft, using hnRNP A/B antibody (Cat No. C15410263) at a 1:100 dilution.

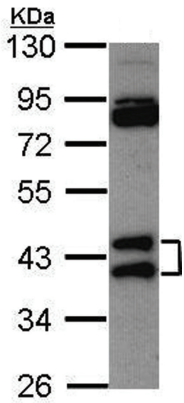
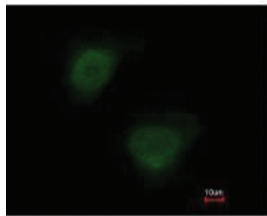


Figure 2. Western blot

Sample: 30 ug of HeLa whole cell lysate
10% SDS PAGE
C15410263 diluted 1:1,000



Costained with Hoechst 33342

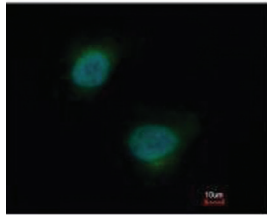


Figure 3. IF

Immunofluorescence analysis of paraformaldehyde-fixed HeLa cells using hnRNP A/B antibody (Cat. No. C15410263) at a 1:200 dilution.

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 12, 2014