

ZNF189 polyclonal antibody - Classic

Cat. No. C15410261

Type: Polyclonal

Source: Rabbit

Lot #: 40730

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	1:100-1:2,000	Figure 2
Western blot	1:500-1:3,000	Figure 1

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

Results

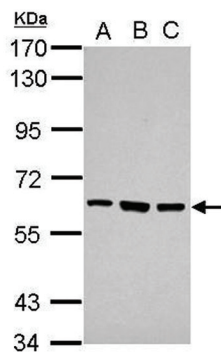


Figure 1. Western Blot

Sample (30 µg of whole cell lysate)

A: Jurkat

B: Raji

C: K562

7.5% SDS PAGE

C15410261 diluted 1:1000

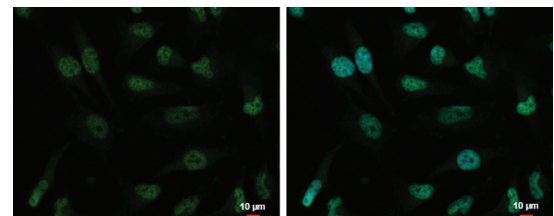
Figure 2. IF

ZNF189 antibody detects ZNF189 protein at nucleus by immunofluorescent analysis.

Sample: HeLa cells were fixed in 4% paraformaldehyde at RT for 15 min.

Green: ZNF189 protein stained by ZNF189 antibody (C15410261) diluted 1:2000.

Blue: Hoechst 33342 staining.



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