

## ASH1L polyclonal antibody - Classic

**Cat. No.** C15410304

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 001

**Size:** 100 µl

**Concentration:** 1 µg/µl

**Specificity:** Human, mouse, rat

**Purity:** Purified by peptide affinity chromatography.

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

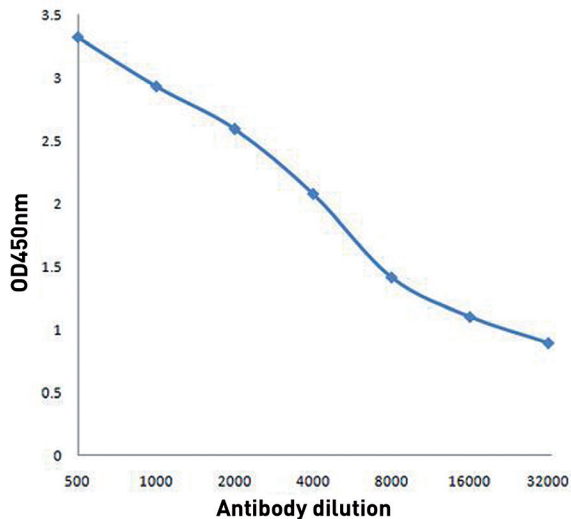
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

### Applications

	Suggested dilution*	Results
ELISA	1:500-1,000	Figure 1
IHC	1:100-500	
IF	1:100-500	

\*Optimal dilution should be determined by the end-user.

### Results



**Figure 1. ELISA results**

ELISA analysis with 2 µg/100 µl antigen and serially diluted ASH1L antibody (1:500, 1:1,000, 1:2,000, 1:4,000, 1:8,000, 1:16,000, 1:32,000).

#### 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 9, 2014