

## ASH2 polyclonal antibody - Pioneer

Cat. No. C15410026

Type: Polyclonal	Specificity: Human
Size: 50 µg	Isotype: NA
Concentration: 2.0 µg/µl	Source: Rabbit
Lot No.: 001	Purity: Protein G purified
Storage buffer: PBS containing 0.05% azide and 0.05% ProClin 300.	Storage conditions: 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.	

Last Data Sheet Update: April 25, 2018

### Description

Alternative names: **ASH2L, ASH2L1, ASH2L2, Bre2**

Polyclonal antibody raised in rabbit against human Ash2 (absent, small, or homeotic 2), using a recombinant protein.

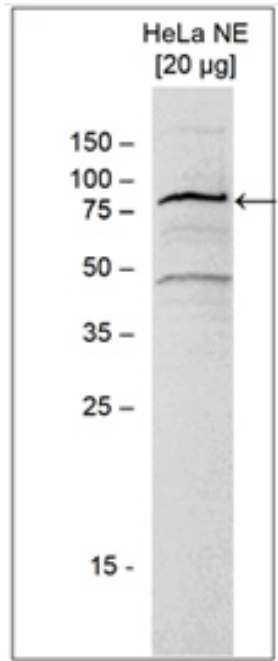
### Applications

Applications	Suggested dilution	References
Western Blotting	1:1,000	Fig 1

### Target Description

ASH2 (UniProtKB/Swiss-Prot entry Q9UBL3) is a component of the SET1/ASH2 histone methyltransferase (HMT) complex. This complex specifically methylates K4 of histone H3, thereby activating transcription. Methylation of K4 is blocked by premethylation of the neighboring K9, a repressor of transcription. This indicates that the SET1/ASH2 HMT complex mediates the crosstalk between K9 methylation and K4 methylation. ASH2 plays a role in hematopoiesis and may be associated with some kinds of leukemia.

**Validation Data**



**Figure 1. Western blot analysis using the Diagenode antibody directed against ASH2**

Western blot was performed on HeLa nuclear extract (HeLa NE, 20 µg) using the Diagenode antibody against ASH2 (Cat. No. C15410026) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest is indicated on the right.