

Anti-acetyl Histone H3 (Lys27), mouse monoclonal antibody

Catalog No. **MA309B**
Lot No. **09001**
Storage **Store below -20 °C**
Expiration Date **Mar 2011**

Product Description

| | |
|---------------|---|
| Clone No. | MABI0309 (CMA309) |
| Host | Mouse |
| Source | Culture supernatant of serum free medium |
| Isotype | IgG1 |
| Size | 100 µl |
| Concentration | 1 mg/ml |
| Antigen | 19 amino acid residues around the Lys27 of human Histone H3.1 |
| Buffer | PBS + 0.05% sodium azaid |

Application notes

| | |
|---------------------|-----------------------|
| Immunoprecipitation | 1-5 µg/5 µl Sepharose |
| Immunostaining | 0.5-1 µg/ml |
| Immunoblot | 0.5-1 µg/ml |

Attention

Mouse IgG has weak combination with protein A and protein G. Please use the beads which connect anti-mouse IgG antibodies such as anti-mouse IgG sepharose for immunoprecipitation.

Reference

Kimura H, Hayashi-Takanaka Y, Goto Y, Takizawa N, Nozaki N. The organization of histone H3 modifications as revealed by a panel of specific monoclonal antibodies. Cell Struct Funct. 2008;33(1):61-73.

MABI 0309

