

## Anti-acetyl Histone H3 (Lys9/27), mouse monoclonal antibody

**Catalog No.**      **MA310B**  
**Lot No.**            **09001**  
**Storage**            **Store below -20 °C**  
**Expiration Date**   **Mar 2011**

### Product Description

|               |  |
|---------------|--|
| Clone No.     | MABI0310 (CMA310)  |
| Host          | Mouse  |
| Source        | Culture supernatant of serum free medium                         |
| Isotype       | IgG2a  |
| Size          | 100 µl   |
| Concentration | 1 mg/ml  |
| Antigen       | 13 amino acid residues from the N-terminus 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 0310

