# Anti-trimethyl Histone H3 (Lys27),

mouse monoclonal antibody

Catalog No. MA323B

Lot No. 11001

Storage Store below -20 °C

Expiration Date Apr 2014

## **Product Description**

Clone No. MABI0323 Host Mouse

Source Culture supernatant of serum free medium

 $\begin{array}{ccc} Isotype & IgG1 \\ Size & 100 \ \mu l \\ Concentration & 1 \ mg/ml \end{array}$ 

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

 $\begin{array}{ll} Immunostaining & 0.5\mbox{-}1\ \mu g/ml \\ Immunoblot & 0.5\mbox{-}1\ \mu g/ml \end{array}$ 

#### 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 0323

