

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

Catalog No. **MA305B**
Lot No. **09001**
Storage **Store below -20 °C**
Expiration Date **Mar 2011**

Product Description

Clone No.	MABI0305 (CMA305)
Host	Mouse
Source	Culture supernatant of serum free medium
Isotype	IgG2b
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 0305

