



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### Anti-H3K27ac monoclonal antibody Catalog #: 25238

<b>Lot #:</b> 141211	<b>Host Species:</b> Mouse
<b>Conc.:</b> 1 µg/µl	<b>Species Reactivity:</b> Human
<b>Size:</b> 50 µg	<b>Immunogen:</b> Synthetic peptide
<b>Clonality:</b> Monoclonal	<b>Purification:</b> Protein A purified

**Description:** Monoclonal antibody raised in Mouse against histone H3 acetylated at lysine 27 (H3K27ac), using a KLH-conjugated synthetic peptide

**Background:**

**Formulation:** PBS containing 0.05% azide and 0.05% ProClin 300

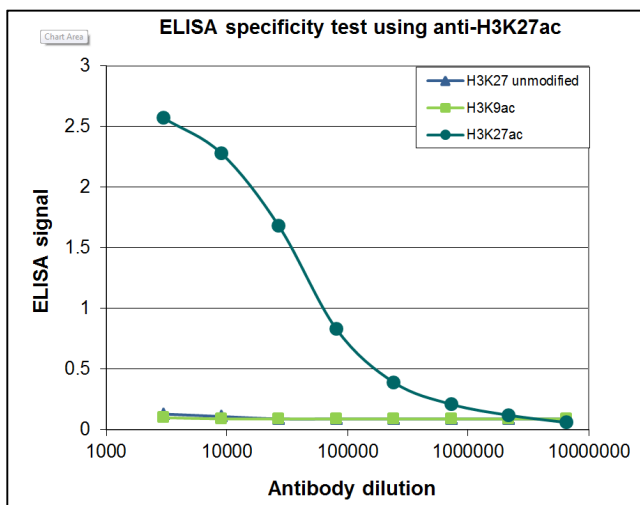
**Applications:** ChIP (1 - 5 µg/ChIP), ELISA (1:3000), WB (1:1000 - 1:2000), IF (1:500)

**Storage/Stability:** Store at -80°C for up to 2 years. Centrifuge after first thaw to maximize product recovery. Aliquot to avoid repeated freeze/thaw cycles. Aliquots may be stored at -20°C for at least one month.

**Warnings:** Avoid freeze/thaw cycles

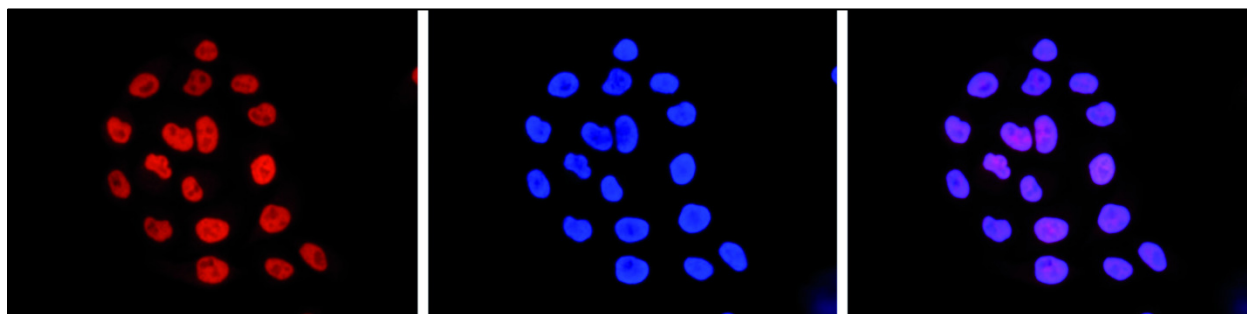
**Notes:** The optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-5 µg per IP

## Quality Assurance:



### Cross reactivity of the monoclonal antibody directed against H3K27ac.

To test the specificity an ELISA was performed using a serial dilution of the monoclonal antibody against H3K27ac (Cat. No. 25238). The wells were coated with peptides containing the unmodified H3K27 region as well as the acetylated H3K27 and the acetylated H3K9. Figure 1 shows a high specificity of the antibody for the peptide containing the modification of interest.



### Immunofluorescence using the monoclonal antibody directed against H3K27ac.

HeLa cells were stained with the antibody against H3K27ac (Cat. No. 25238) and with DAPI. Cells were fixed with 4% formaldehyde for 10 minutes and blocked with PBS/TX-100 containing 5% normal goat serum and 1% BSA. The cells were immunofluorescently labelled with the H3K27ac antibody (left) diluted 1:500 in blocking solution followed by an anti-mouse antibody conjugated to Alexa594. The middle panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.