



6405 Mira Mesa Blvd, Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data Sheet

Anti-H3K27ac monoclonal antibody

Catalog #: 25238

Lot #: 220324	Host Species: Mouse
Conc.: 1 µg/µl	Species Reactivity: Human
Size: 50 µg	Immunogen: Synthetic peptide
Clonality: Monoclonal	Purification: Protein A purified

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

Formulation: PBS containing 0.05% azide

Applications: ChIP (1 - 2 µg per IP), ELISA (1:3000), WB (1:1000 - 1:2000), IF (1:500)

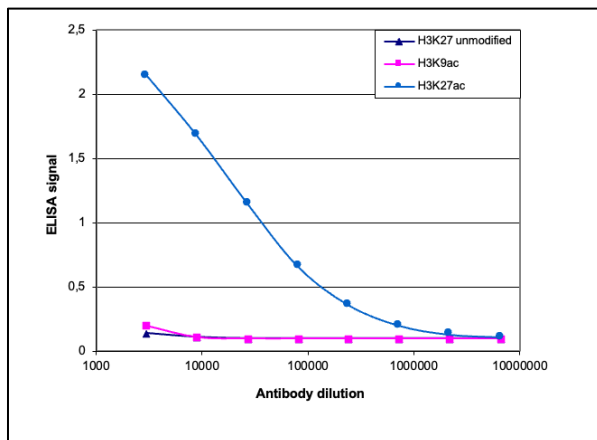
Storage/Stability: Store at -80°C for up to 2 years. Centrifuge after first thaw to maximize product recovery. Aliquot to avoid repeated freeze/thaw cycles. Aliquots may be stored at -20°C for at least one month.

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

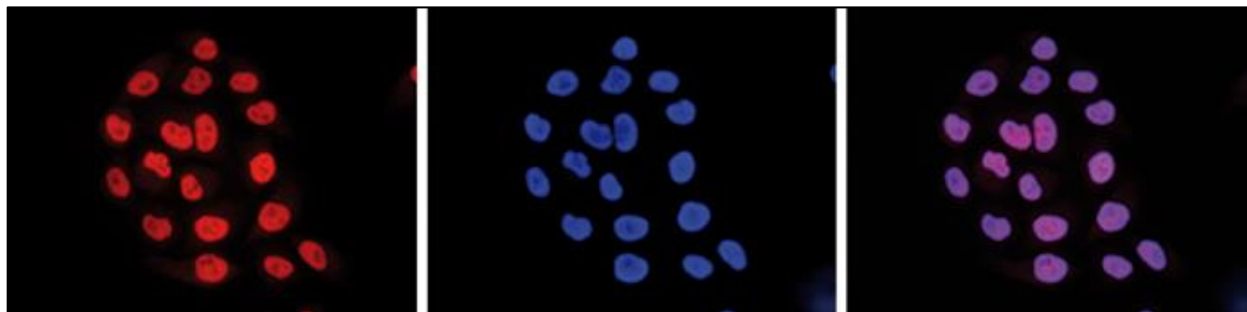
Quality Assurance:

ELISA specificity test using monoclonal antibody against H3K27ac



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.