



Histone H4 acetylated Monoclonal Antibody

Cat No: G01058 - 100 μ L

General Data

Shipping:	wet ice
Stability:	Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative
Formulation:	Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles
Host:	Mouse
Antigen:	synthetic peptide: MSGRG(acetyl-K)GG(acetyl-K)GLG(acetyl-K)GGA(acetyl-K)RHRKVL(R)-C corresponding to aa 1 to 24 of human histone H4. This sequence is conserved in a wide range of species.
Clone:	2C2
Isotype:	IgG1, κ
Application(s):	ELISA Western Blot Immunoprecipitation Immunofluorescence Recommended dilutions: 1/500-1/5000
Specificity:	Acetylated histone H4 at positions K5, K8, K12 or K16



Product Overview

Histone H4 is a structural component of the nucleosome and is subject to covalent modification including acetylation and methylation, which may alter expression of genes located on DNA associated with its parent histone octamer. The Histone H4 clone 3HH-2C2 monoclonal recognize H4 when actelylated on at list on of the for Lysine residues K5, K8, K12 or K16.

FP/20/20

For research laboratory use only – Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux - France
Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>
<https://www.bertin-bioreagent.com>



Histone H4 acetylated Monoclonal Antibody

Cat No: G01058 - 100 ÅµL

Scientific Literature

Brand M, Moggs JG, Oulad-Abdelghani M, Lejeune F, Dilworth FJ, et al. (2001) UVdamaged DNA-binding protein in the TFTC complex links DNA damage recognition to nucleosome acetylation. *Embo J* 20: 3187-3196.

Dilworth FJ, Fromental-Romain C, Yamamoto K, Chamb

FP/20/20

For research laboratory use only – Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux - France
Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>
<https://www.bertin-bioreagent.com>