

Specifications (926-42218)

- Size: 100 µL
- Storage: -20 °C
- Species Cross-Reactivity: human, mouse, rat, monkey
- Target Molecular Weight: 17 kDa
- Isotype: Mouse IgG₃
- Specificity/Sensitivity: Detects endogenous levels of total Histone H3 protein.
- Immunogen: A peptide that is specific to the carboxy terminus of human histone H3 protein
- Storage Buffer: 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg/mL BSA, 50% glycerol, and <0.02% sodium azide

Note: Do not aliquot the antibody.

Warning: Sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.

- Tested Application: Western blot, Immunohistochemistry, Immunofluorescence, Flow Cytometry, Chromatin Immunoprecipitation

Recommended Dilution

- Western Blotting: 1:1000

Applications

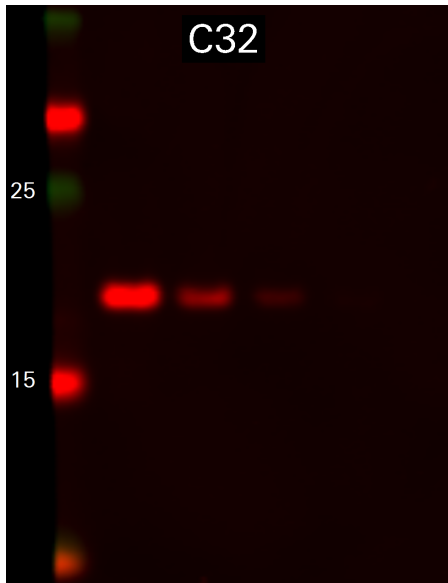
The Histone H3 primary antibody can be used as an internal loading control for normalization and is particularly effective when detecting target proteins in nuclear extracts.

The expression of Histone H3, or any housekeeping protein (HKP), should be validated to ensure that its expression does not change under experimental conditions.

Once validated, Histone H3 primary antibodies can be used for the detection of Histone H3 when performing multiplex Western blot detection.

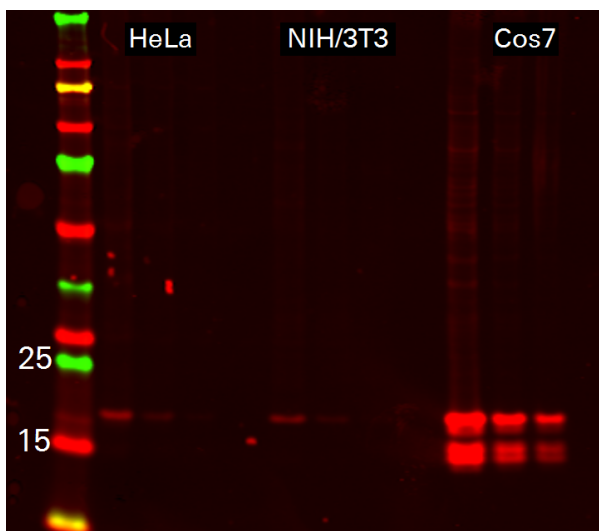
Detect Histone H3 Mouse Monoclonal Antibody with IRDye® Goat anti-Mouse or IRDye Donkey anti-Mouse secondary antibodies.

Histone H3 Mouse Monoclonal Antibody in C32 Lysates



Histone H3 Mouse Monoclonal Antibody was detected in C32 cells. C32 lysates were diluted from 2.5 μ g to 156 ng. Lysates were separated on 4-12% Bis-Tris Gels electrophoresed at 200V for 45 minutes in MES Running Buffer and transferred to nitrocellulose membranes in Tris Glycine buffer at 100V for 65 minutes. Blots were blocked in Intercept (PBS) Blocking Buffer and probed with Histone H3 Mouse Monoclonal Antibody and detected on an Odyssey CLx Imager.

Histone H3 Mouse Monoclonal Antibody in HeLa, NIH/3T3, and Cos7 Lysates



Histone H3 Mouse Monoclonal Antibody detected in HeLa, NIH/3T3, and COS7 lysates. Lysates were diluted from 2.5 μ g to 156 ng. Lysates were separated on 4-12% Bis-Tris Gels electrophoresed at 200V for 45 minutes in MES Running Buffer and transferred to nitrocellulose membranes in Tris Glycine buffer at 100V for 65 minutes. Blots were blocked in Intercept (PBS) Protein-Free Blocking Buffer and probed with Histone H3 Mouse Monoclonal Antibody and detected on an Odyssey CLx Imager.

Limitation of Liability and Limited Use Label License

For Research Use Only. Not intended for human therapeutic or diagnostic use.

By using this product, you agree to the Limitation of Liability and the Limited Use Label License available online at licor.com/packinserts under the **Pack Inserts** heading (the "License"). The License is incorporated herein by this reference, and the License may be updated from time to time.



© 2020 LI-COR, Inc. LI-COR, In-Cell Western, Chameleon, Intercept, IRDye, and Odyssey are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com