

## Specifications (926-42216)

- Size: 100  $\mu$ L
- Storage: -20 °C
- Species Cross-Reactivity: human, mouse, rat, monkey
- Target Molecular Weight: 37 kDa
- Isotype: Rabbit IgG
- Specificity/Sensitivity: Detects endogenous levels of total GAPDH protein. May cross-react with pig.
- Immunogen: A synthetic peptide that corresponds to residues near the carboxy terminus of human GAPDH
- Storage Buffer: 10 mM HEPES (pH 7.5), 150 mM NaCl, 100  $\mu$ g/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 Applications: Western blot, Immunohistochemistry, Immunofluorescence

## Recommended Dilution

- Western Blotting: 1:1000

## Applications

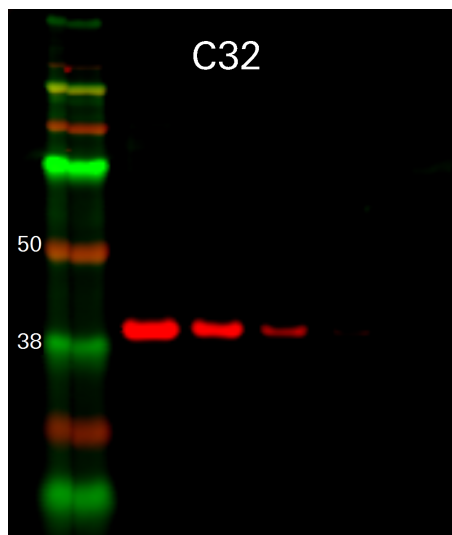
Glyceraldehyde-3-phosphate dehydrogenase (GAPDH) is a constitutively expressed housekeeping protein (HKP). The GAPDH primary antibody can be used as an internal loading control for normalization.

The expression of GAPDH, or any HKP, should be validated to ensure that its expression does not change under experimental conditions.

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

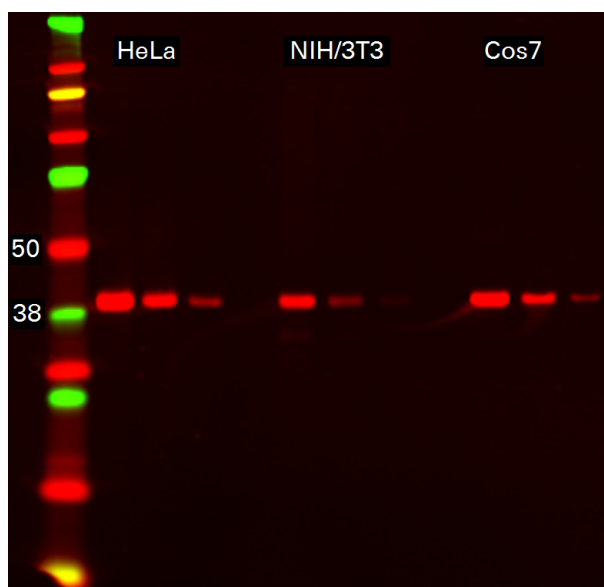
Detect GAPDH Rabbit Monoclonal Antibody with IRDye® Goat anti-Rabbit or IRDye Donkey anti-Rabbit secondary antibodies.

## GAPDH Rabbit Monoclonal Antibody in C32 Lysates



**GAPDH Rabbit Monoclonal Antibody was detected in C32 cells.** C32 lysates were diluted from 2.5  $\mu$ 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 GAPDH Rabbit Monoclonal Antibody and detected on an Odyssey CLx Imager.

## GAPDH Rabbit Monoclonal Antibody in HeLa, NIH/3T3, and Cos7 Lysates



**GAPDH Rabbit Monoclonal Antibody detected in HeLa, NIH/3T3, and COS7 lysates.** Lysates were diluted from 2.5  $\mu$ 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 GAPDH Rabbit 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](http://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](mailto:biosales@licor.com)  
[licor.com/bio](http://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](mailto: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](mailto:bio-eu@licor.com)