

Mouse Anti- β -Actin Monoclonal Antibody

CATALOG NUMBER:	AKR-002	STORAGE:	-20°C
QUANTITY AND CONCENTRATION:	100 μ g of Protein A purified antibody at 1 mg/mL in 10 mM PBS, pH 7.2, containing 0.05% NaN ₃		
SHELF LIFE:	1 year from date of receipt under proper storage conditions; aliquot to avoid multiple freeze thaw cycles		
HOST SPECIES:	Mouse		
ISOTYPE:	IgG2b		
CLONE:	BA3R		
IMMUNOGEN:	N-terminal β -actin conjugated to KLH		
SPECIFICITY:	Native and denatured β -actin (~42 kD)		
REACTIVITY:	Human, mouse, rat, rabbit, chicken; other species not tested		
APPLICATION:	Immunoblot (1:1000 to 1:10,000; incubate 1 hr at room temp) Immunostaining (1:500 to 1:2000) ELISA or Dot blot (user to determine working dilution)		

Example of Results

The following figures demonstrate typical results. One should use the data below for reference only. This data should not be used to interpret actual results.

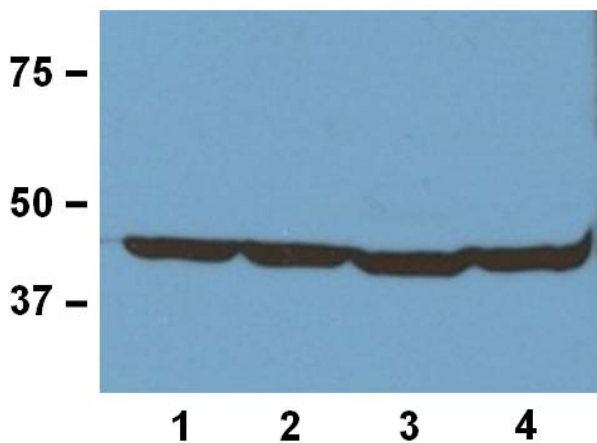


Figure 1. Immunoblotting with Anti- β -Actin clone #BA3R. Anti- β -Actin antibody at 1 μ g/mL (1:1000 dilution) at 20 μ g/lane tissue lysate from human (1), mouse (2), rat (3), and rabbit (4).

Recent Product Citation

Orlandi, A. et al. (2015). ERCC1 uinduction after oxaliplatin exposure may depend on KRAS mutational status in colorectal cancer cell line: in vitro veritas. *J Cancer*. **6**:70-81.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2011-2015: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.