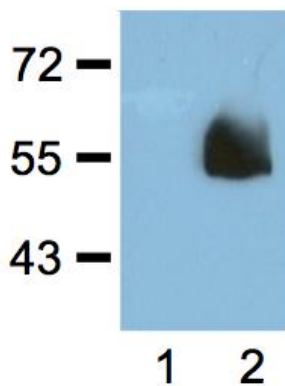


## Mouse Anti-HA Tag Monoclonal Antibody

<b>CATALOG NUMBER:</b>	AKR-006	<b>STORAGE:</b>	-20°C
<b>QUANTITY AND CONCENTRATION:</b>	100 µg of Protein A purified antibody at 1 mg/mL in 10 mM PBS, pH 7.2, containing 0.05% NaN <sub>3</sub>		
<b>SHELF LIFE:</b>	1 year from date of receipt under proper storage conditions; aliquot to avoid multiple freeze thaw cycles		
<b>HOST SPECIES:</b>	Mouse		
<b>ISOTYPE:</b>	IgG3		
<b>CLONE:</b>	HA.C5		
<b>IMMUNOGEN:</b>	YPYDVDPDYA synthetic peptide conjugated to KLH		
<b>SPECIFICITY:</b>	Recognizes the epitope of N-terminal, C-terminal, or internal YPYDVDPDYA tags		
<b>SENSITIVITY:</b>	1-5 ng of HA-tagged fusion proteins (immunoblot/ECL)		
<b>APPLICATION:</b>	Immunoblot (1:1000 to 1:3000; incubate 1 hr at room temp) Immunostaining (1:500 to 1:2000) ELISA, Dot blot or Immunoprecipitation (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-HA Tag clone #HA.C5.** Anti-HA Tag antibody at 1 µg/mL (1:1000 dilution). **Lane 1:** Untransfected control. **Lane 2:** HEK293 cells transfected with HA-tagged protein vector.

## **Recent Product Citations**

1. Mohammed, S. et al. (2020). Regulatory role of SphK1 in TLR7/9-dependent type I interferon response and autoimmunity. *FASEB J.* doi: 10.1096/fj.201902847R.
2. Starostová, M. et al. (2014). The oncoprotein v-Myb activates transcription of Gremlin 2 during in vitro differentiation of the chicken neural crest to melanoblasts. *Gene.* **540**:122-129.

## **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](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2011-2020: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.