

Anti-HACE1 Recombinant Mouse Monoclonal Antibody

Store at 2°C to 8°C (short-term), or -20°C (long-term)

Catalog Number 730103

Pub. No. MAN0007644 Rev. 1.01

Clonality: Monoclonal	Quantity: 100 µg	Volume: 200 µL	Concentration: 0.5 mg/mL
Host/Class: Mouse IgG	Reactivity: Human HACE1	Predicted Reactivity: Human	

Product Description

HACE1 is an E3 ubiquitin-protein ligase involved in Golgi membrane fusion and regulation of small GTPases. It is recruited to Golgi membranes through interaction with Rab proteins, and acts as a regulator of Golgi membrane formation by mediating ubiquitination during mitotic Golgi disassembly. Ubiquitination acts as a regulatory signal for postmitotic Golgi membrane fusion. HACE1 also plays a role in host defense by mediating the ubiquitination of active GTP-bound RAC1 to catalyze its degradation. RAC1 degradation is important during pathogen defense, as RAC1 is targeted by a number of bacterial toxins. HACE1 may also act in transcription regulation through interaction with RAR beta.

This antibody was developed against a fully folded protein domain, so it does not show good reactivity against denatured proteins, and may not work in western blot applications.

Product Specifications

Immunogen:	Folded protein domain corresponding to amino acids 420–909 (HECT domain) of human HACE1.
Apparent MW:	102.3 kDa
Gene ID:	57531
Protein Accession No.:	Q81YU2
Sequence Identity:	Human
Isotype:	IgG ₁
Lot:	See product label

Product Applications

Application	Species	Test Material	Concentration
Immunoprecipitation + Mass Spectrometry	Human	HEK293 cells	2 µg per IP

Storage and handling

Store the antibody at 2°C to 8°C for up to 1 month, or -20°C for long storage. Avoid repeated freezing and thawing.

Stability

When stored as instructed, expires one year from date of receipt unless otherwise indicated on product label.

Storage Buffer

Phosphate buffered saline (PBS) with 0.09% sodium azide.



CAUTION! Sodium azide is extremely toxic and may react with lead and copper plumbing to form highly explosive metal azides. Properly dispose of solutions containing sodium azide. Read the Safety Data Sheet (SDS) and follow the handling instructions. Wear appropriate protective eyewear, clothing and gloves. SDSs are available at www.lifetechnologies.com/support.

Product Documentation

To obtain a Certificate of Analysis or Safety Data Sheet (SDS), visit <http://www.lifetechnologies.com/support>.








Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Related Products

Product Name	Quantity	Catalog No.
iBlot [®] Dry Blotting System	1 unit	IB1001
WesternBreeze [®] Chromogenic Kit Anti-Rabbit	1 kit	WB7105
WesternBreeze [®] Chemiluminescent Kit, Anti-Rabbit	1 kit	WB7106
Goat anti-mouse (H+L), HRP conj.	1 mg	G21040
Goat anti-rabbit (H+L), HRP conj.	1 mg	G21234
Goat anti-mouse (H+L), AP conj.	1 mg	G21060
Goat anti-rabbit (H+L), AP conj.	1 mg	G21079
Streptavidin Agarose	5 mL	SA10004

Explanation of Symbols

Symbol	Description	Symbol	Description	Symbol	Description
	Manufacturer		Catalog number		Batch code
	Use by		Temperature limitation		
	Consult instructions for use		Caution, consult accompanying documents		

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

NOTICE TO PURCHASER: Limited Use Label License No. 327: Recombinant Antibody Technology: The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research and manufacturing conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product, (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components for contract manufacturing services; (2) use of the product or its components to provide a service, information, or data; (3) use of the product itself or its components as a therapeutic, diagnostic or prophylactic; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research or manufacturing. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies is willing to accept return of the product for a full refund. For information on obtaining additional rights, please contact outlicensing@lifetech.com.

© 2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com