

BRF1/2 Antibody

Publication Part no. MAN0004925

Rev. Date: 21 July 2011

Catalog Number: A13961

Store at -20°C

Clonality: Polyclonal
Volume: 100 µL

Host/Class: Rabbit IgG
Reactivity: Human BRF1/2
Predicted Reactivity: Human, Mouse, Monkey, Rat

Product Description

BRF1 (butyrate response factor 1) and BRF2 (butyrate response factor 2) belong to the TIS11 family of CCCH zinc-finger proteins. This family of proteins bind to AU-rich elements (ARE) found in the 3'-untranslated regions of mRNAs and promote de-adenylation and rapid degradation by the exosome. These proteins play a critical role in cell growth control by regulating the mRNA turnover of multiple cytokines, growth factors and cell cycle regulators, including GM-CSF, TNF α , IL-2, IL-3 and IL-6. Insulin-induced stabilization of ARE-containing transcripts is mediated by Akt/PKB phosphorylation of BRF1 at Ser⁹², which results in binding by 14-3-3 protein and inactivation of BRF1.

Product Specifications

Immunogen: Synthetic peptide corresponding to the carboxy terminus of human BRF1 protein

Alternate Names: BRF, ERF1/2, EGF response factor 1/2, GTF3B, TAF3B2, TAF3C, TAFIII90, TF3B90, TFIIB90, TIS11B/D, hBRF, ZFP36L1/2

Apparent MW: 40-50, 62 kDa

Gene ID: 677

Protein Accession No.: Q07352

Sequence Identity: Human

Lot: See product label

Product Applications

Application	Species	Test Material	Concentration
Western blotting	Human	HCT 15 cells	1:1000
Immuno fluorescence (IF-IC)	Human	DLD-1 cells	1:25
Flow Cytometry	Human	A549 cells	1:100

Storage and Handling

Store the antibody at -20°C. Do not make aliquots of the antibody.

Stability

Expires one year from date of receipt when stored as instructed.

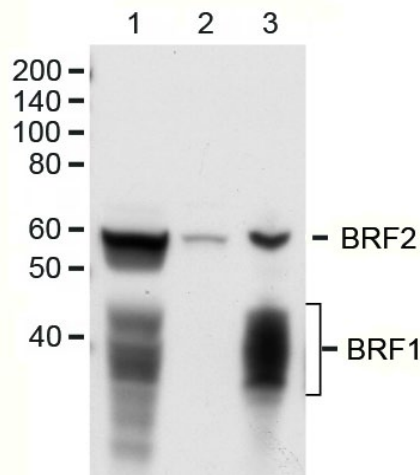


Figure 1 Western blot analysis of cell lysates from HT1080 (lane 1), HT1080 BRF1 -/- (lane 2), and HCT15 (lane 3) cells, using BRF1/2 Antibody.

For research use only. Not for human or animal therapeutic or diagnostic use.

Manufacturing Site • 7335 Executive Way • Frederick • MD 21704 • E-mail: techsupport@invitrogen.com

Storage Buffer

10 mM sodium HEPES (pH 7.5), 150 mM NaCl, 100 µg/mL BSA and 50% glycerol.

Safety Data Sheets (SDS)

Safety Data Sheets (SDSs) are available at www.invitrogen.com/sds.









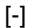
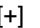


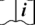
Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.invitrogen.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

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
Nitrocellulose, 0.2 µm	20/pack	LC2000

Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		<i>In vitro</i> diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.invitrogen.com/support or email techsupport@invitrogen.com

www.lifetechnologies.com

 life
technologies™