

## GABA Transporter (GAT) 1, Rabbit Polyclonal Antibody

Catalog Number: P21948

Store at -20°C

**Clonality:** Polyclonal  
**Volume:** 100 µL

**Host/Class:** Rabbit  
**Reactivity:** GABA transporter 1  
**Predicted Reactivity:** Rat

### Product Description

GABA plasma membrane transporters (GATs) influence synaptic neurotransmission by high-affinity uptake and release of Gamma-aminobutyric acid (GABA). Four distinct GABA transporters have been identified to date: GAT-1, GAT-2, GAT-3, and BGT-1. GAT-1, the most abundant of the transporters, is found predominantly in neurons, but also in some specialized glia. GAT-1 is thought to play a key role in epileptogenesis.

### Product Specifications

**Immunogen:** Peptide corresponding to amino acid residues from the C-terminal region of rat GAT-1.  
**Apparent MW:** ~67 kDa  
**Gene ID:** 79212  
**Protein Accession No.:** P23978  
**Sequence Identity:** Rat  
**Sequence Homology:** Mouse  
**Lot:** See product label

### Product Applications

Application	Dilution
Western blotting	1:1000
Immunohistochemistry	1:100-1:200

### Storage and Handling

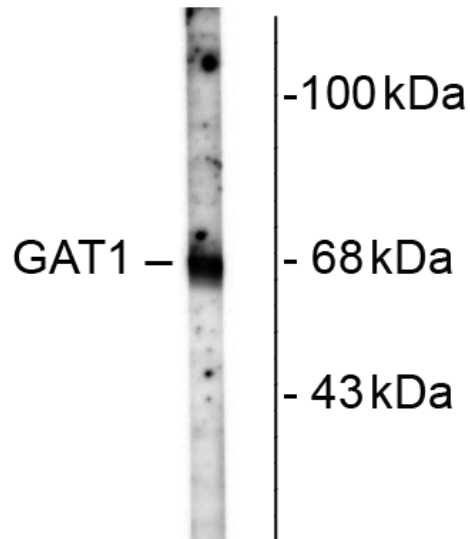
Store the antibody at -20°C. **Avoid freeze/thaw cycles.**

### Stability

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

### Storage Buffer

100 µL in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg per mL BSA and 50% glycerol.



**Figure 1** Western blot of rat hippocampal homogenate showing specific immunolabeling of the ~67 kDa GAT-1 protein.

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

E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

Life Technologies Corporate Headquarters • 5791 Van Allen Way • Carlsbad • CA 92008 • Tel: 800.955.6288 • [www.invitrogen.com](http://www.invitrogen.com)

MAN0003834

Rev. Date: 21 March 2011

## Safety Data Sheets (SDS)

Safety Data Sheets (SDSs) are available at [www.invitrogen.com/sds](http://www.invitrogen.com/sds).








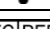

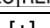
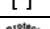
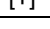

## Certificate of Analysis

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](http://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](mailto: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.