

## Datasheet for

**Catalog number:**

**Notice:** Undefined index: quantity in  
[/home/genec2/httpdocs/product/antibody/antibody\\_pdf.php](/home/genec2/httpdocs/product/antibody/antibody_pdf.php) on line **44**

**Background:**

■ **Product Information**

**Symbol:**

**Alias:**

**Cross reactivity:**

**Expected Protein Weight:**

**Speices:**

**Source:**

**Isotype:**

**Purity:**

**Storage:**

■ **Immunogen Information**

**Gene ID:**

**Swiss:**

**Immunogen:**

■ **Recommended Dilutions**

**Notice:** Undefined offset: 1 in  
[/home/genec2/httpdocs/product/antibody/antibody\\_pdf.php](/home/genec2/httpdocs/product/antibody/antibody_pdf.php)  
on line **132**

**Citation of product:** If use of this item results in a publication, please use this information: (

**Notice:** Undefined index: quantity in

[/home/genec2/httpdocs/product/antibody/antibody\\_pdf.php](/home/genec2/httpdocs/product/antibody/antibody_pdf.php) on line **138**

-; Genecopoeia, Inc, Rockville, MD).

GeneCopoeia Products are for Research Use Only  
9620 Medical Center Drive, #101 Rockville, MD 20850 USA  
Tel: 301-762-0888:Fax:301-762-3888  
Email: [support@genecopoeia.com](mailto:support@genecopoeia.com); Web:[www.genecopoeia.com](http://www.genecopoeia.com)

Copyright © 2024 GeneCopoeia, Inc.