

Datasheet for

Catalog number:

Notice: Undefined index: quantity in
/home/genec2/httpdocs/product/antibody/protein_pdf.php on line
54
-

Description:

■ **Protein Information**

Species:

AA sequence:

Uniprot ID:

Expression host:

Tag:

Calculated MW:

Protein families:

Protein pathways:

■ **Reference Information**

Symbol:

Alias:

Accession:

Gene ID:

RefSeq:

RefSeq size:

RefSeq ORF:

Cytogenetics:

■ **Product Information**

Concentration:

Purity

Buffer:

Shipping:

Storage:

Stability:

Citation of product: If use of this item results in a publication, please use this information: (; 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; Web: www.genecopoeia.com

Copyright © 2024 GeneCopoeia, Inc.