

## Datasheet for HRP conjugated Rabbit Anti-Monkey IgD mAb

**Catalog number:** mAb-02619-100

**Background:** Secondary antibodies are affinity-purified antibodies which will work with target-specific primary antibody in the detection, sorting or purification of its specified target. Secondary antibodies offer increased versatility enabling users to use many detection systems (e.g. HRP, AP, fluorescence). They can also provide greater sensitivity through signal amplification as multiple secondary antibodies. Most commonly, secondary antibodies are generated by immunizing the host animal (different from host species of primary antibody) with a pooled population of normal immunoglobulins from the host species of primary antibody and can be further purified and modified (i.e. antibody fragmentation, label conjugation, etc.) to ensure well-characterized specificity to corresponding normal immunoglobulins.

### ■ Product Information

**Symbol:** Monkey IgD  
**Alias:**  
**Cross reactivity:** Cynomolgus monkey  
**Expected Protein Weight:**  
**Speices:**  
**Source:** Rabbit  
**Isotype:** IgD  
**Purity:** Affinity purification  
**Storage:** Store at -20°C. Avoid freeze / thaw cycles. Buffer: PBS with 0.05% proclin300, 0.05% BSA, 50% glycerol, pH7.3.

### ■ Immunogen Information

**Gene ID:**  
**Swiss:**  
**Immunogen:** Monkey IgD

### ■ Recommended Dilutions

**ELISA** 1:5000 -1:10000

**Citation of product:** If use of this item results in a publication, please use this information: HRP conjugated Rabbit Anti-Monkey IgD mAb(mAb-02619-100; 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.