



**GeneCopoeia Inc.**

9620 Medical Center Drive, Suite 101

Rockville, MD 20850

USA

Phone: 301-762-0888

Toll free: 1-301-762-0888

Fax: 301-762-3888

Web: [www.genecopoeia.com](http://www.genecopoeia.com)

Inquiry: [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)

Technical Support: [support@genecopoeia.com](mailto:support@genecopoeia.com)