

VitroView™ Histology Services



GeneCopoeia™
Expressway to Discovery

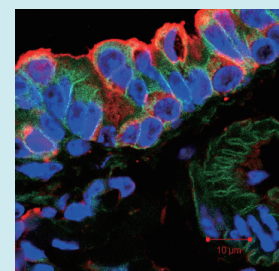
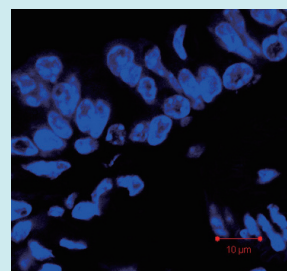
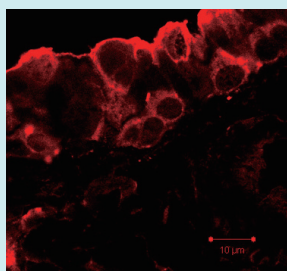
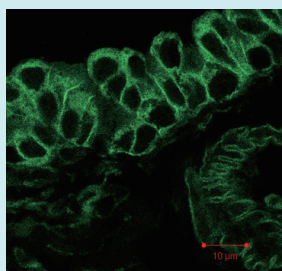
GeneCopoeia provides full histology services with high quality, competitive pricing, and fast turnaround.

Located in Rockville, Maryland, USA, we receive mail orders and also provide FREE PICKUP and DELIVERY for local customers.

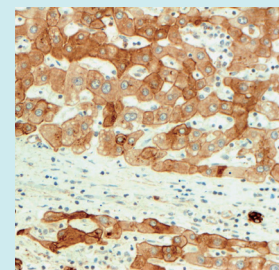
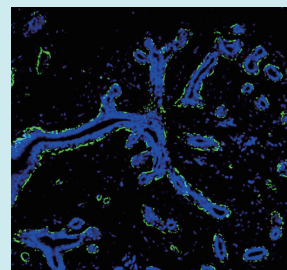
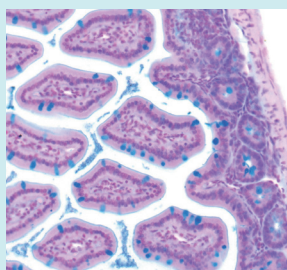
Please visit our website to see more images and details.

FREE PICKUP and FREE DELIVERY
in Bethesda, Rockville, Gaithersburg and Germantown, Maryland

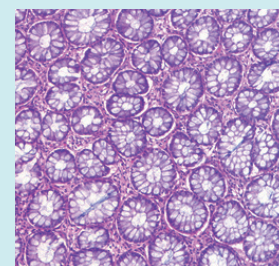
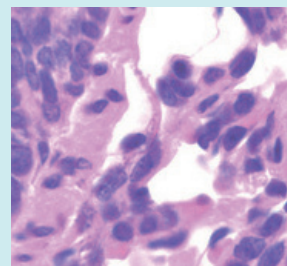
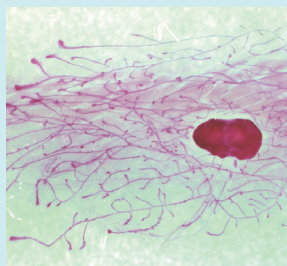
Double IF Staining
(all four)



Alcian Blue Stain *(left)*
Smooth Muscle Actin IF *(middle)*
Pan-keratin IHC/DAB *(right)*



Whole Mount Staining *(left)*
H&E Stain *(middle and right)*



VitroView™ Histology Services

GeneCopoeia provides the following histology services

Routine histology	<ul style="list-style-type: none"> • Whole histological preparation of mammalian species • Paraffin processing • Serial sections • H&E stain
Specialized histology	<ul style="list-style-type: none"> • RNase cutting precaution • Decalcified bone techniques • Special stain
Paraffin-embedding/Sectioning for culture cells	
Cryotomy (frozen sections)	
Immunohistochemistry	
Immunofluorescent (IF) stain (Single or double antibody staining)	
<i>In situ</i> apoptosis (TUNNEL) assay	
<i>In Situ</i> BrdU cell proliferation assay	
Formalin-fixed, paraffin-embedded (FFPE) tissue analysis services	<ul style="list-style-type: none"> • PCR from FFPE tissue/slides • RT-PCR from FFPE tissue/slides • Western blot from FFPE tissue/slides • DNA isolation from FFPE tissue samples • RNA isolation from FFPE tissue samples • Protein extraction from FFPE tissue samples
Mouse genotyping services	<ul style="list-style-type: none"> • Multiple sample types are accepted, including tails, ears or formalin-fixed, paraffin-embedded (FFPE) tissues.

GeneCopoeia, Inc.

9620 Medical Center Drive, Suite 101
Rockville, MD 20850, USA

Email inquiry@genecopoeia.com
Tel +1 (301) 762-0888
Toll free +1 (866) 360-9531
Fax +1 (301) 762-3888
Website www.genecopoeia.com

GeneCopoeia™
Expressway to Discovery

