



## MaxArray™ Human Normal Tissue Microarray Slides

**Qty:** 5 slides

**Catalog No.** 75-4013

**Lot No.**

Invitrogen's MaxArray Human Normal Tissue Arrays contain 30 human tissue samples arrayed on Superfrost Plus Microscope slides. Every package contains 5 unstained slides. Each slide contains 30 different tissues from 30 different organs. Each tissue sample (tissue core) is ~1.5 mm in diameter. The type of tissue and their respective position on the array is indicated in the tissue reference chart. MaxArray tissues are fixed in 10% neutral buffered formalin and embedded in paraffin. Each slide contains a 4 µm tissue section. MaxArray Human Normal Tissue Microarray slides are for research use only.

### STORAGE

Store at room temperature in a cool dry place.

### QUALITY CONTROL

1. Tissues are surgically procured by qualified individuals.
2. H & E slides of donor blocks are examined by Histologists to ensure that the cores are the best representation of subject tissue.

### DEPARAFFINIZATION/REHYDRATION PROTOCOL

1. Deparaffinize slides in xylene for 5 minutes.
2. Place in xylene a second time for an additional 5 minutes.
3. Place the slide in absolute alcohol 2 times for 5 minutes each time.
4. 95% alcohol for 5 minutes.
5. 80% alcohol for 5 minutes.
6. Place slide into distilled water rinse until ready to use.

### IMMUNOHISTOCHEMISTRY (IHC)

1. Once deparaffinized and rehydrated, arrays can be analyzed by traditional IHC methods. Any antibody that has been proven effective in formalin fixed immunohistochemistry or immunocytochemistry should be applicable to MaxArrays.
2. Blocking and incubations conditions are performed with the normal procedure according to the appropriate antibody specifications.
3. If humidity is required, place the slide on a damp paper towel in a plastic container.
4. All wash solutions and reagents should be gently applied to the array slide to prevent dislodging of the cores.

### APPLICATION

MaxArray Human Normal Tissue Microarray slides can be used in many applications including:

1. Immunohistochemistry
2. In situ hybridization, RNA or DNA.
3. Fluorescent in situ hybridization (FISH)\*\* or Chromogenic in situ hybridization (CISH™).
4. MaxArray Human Normal Tissue Microarray slides are for Research Use Only.

\*\*If using fluorescence, please note that tissues inherently autofluoresce and the appropriate filter should be used to distinguish between true fluorescence and autofluorescence. For more information please contact Invitrogen technical support.

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

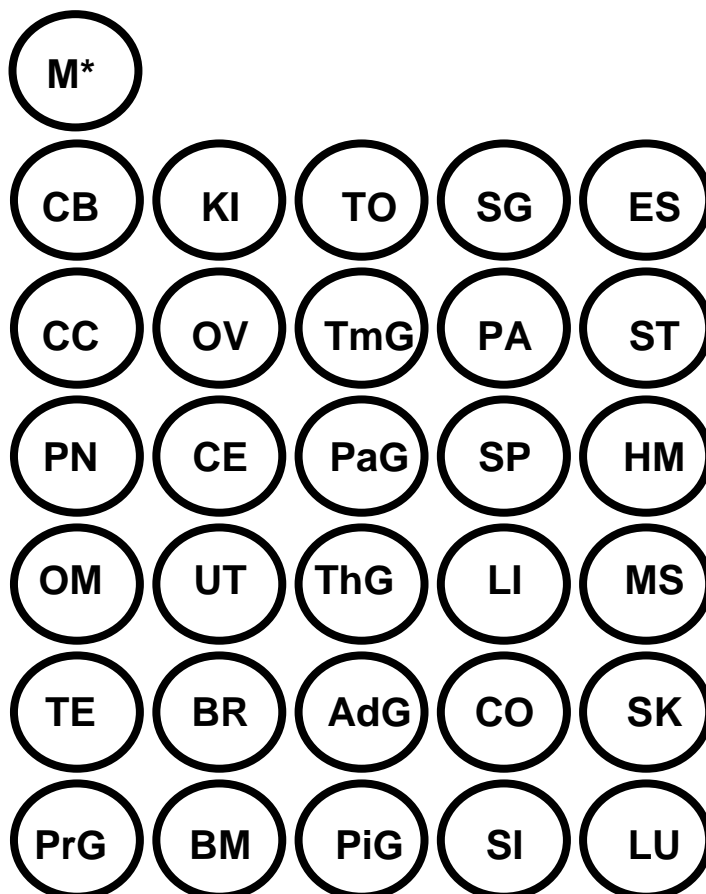
PIN 30970

(Rev 08/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

## TISSUE REFERENCE CHART

Below is a list of the tissues that are included in this array. The respective tissue positions are arranged on the chart in the same position as on the array. Orient yourself by locating the **M\*** (marking tissue) on the array. The marking tissue used is placenta.



### Human Normal Tissue:

LU= Lung  
SK= Skin  
MS= Muscle, skeletal  
HM= Heart muscle  
ST= Stomach  
ES= Esophagus  
SI = Small Intestine  
CO= Colon  
LI= Liver  
SP= Spleen  
PA= Pancreas  
SG= Salivary Gland  
PiG= Pituitary Gland  
AdG= Adrenal Gland  
ThG= Thyroid Gland  
PaG= Parathyroid Gland  
TmG= Thymus Gland  
TO= Tonsil  
BM= Bone marrow  
BR= Breast  
UT= Uterus  
CE= Cervix  
OV= Ovary  
KI= Kidney  
PrG= Prostate Gland  
TE= Testis  
OM= Omentum  
PN= Peripheral Nerve  
CC= Cerebral Cortex  
CB= Cerebellum

**M\*:** Marking tissue (placenta tissue) for orientation of the MaxArray.

MaxArray™, CISH™ and Zymed® are trademarks of Zymed Laboratories, Inc.

**For Research Use Only**

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PIN 30970

(Rev 08/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.