



MaxArray™ Multi-Species (Mouse,Rabbit,Rat,Sheep) Microarray Tissue Slides

Qty: 4 slides + 1 H & E slides

Catalog No. 75-2053

Lot No.

Invitrogen's MaxArray™ Multi-Species Tissue Microarrays contain 60 tissue samples arrayed on Superfrost Plus Microscope slides. Every package contains 4 unstained slides and 1 hematoxylin and eosin (H & E) stained slide for your reference. Each slide contains 15 different tissues from 4 different species (Mouse, Rabbit, Rat and Sheep) for a total of 60 tissue samples per slide. Each tissue sample (tissue core) is 1.5 mm in diameter. The type of tissue and their respective position on the array is indicated below in the tissue reference chart. MaxArray™ tissues are fixed in 10% neutral buffered formalin for 12-24 hrs. and embedded in paraffin. Each slide contains a 4 µm tissue section. MaxArray™ Multi-Species Tissue slides are for research use only.

STORAGE

Store at 2 – 8°C*.

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. Place slide 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 incubation conditons 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 animal tissue 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™ Multi-species tissue 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.

* These recommendations can be only regarded as general guidelines, since different epitopes are known to show diverging susceptibility to duration and conditions of storage.

www.invitrogen.com

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

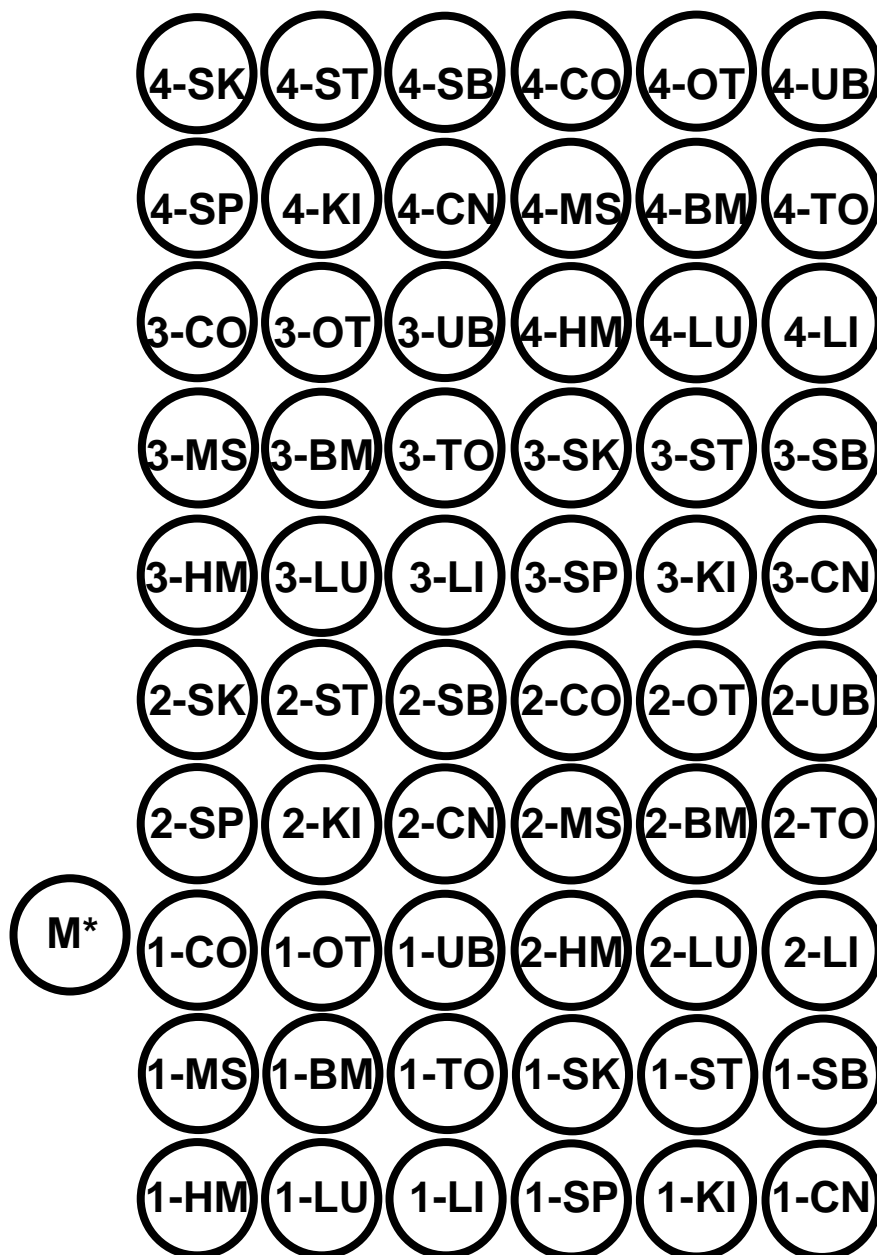
PI 75-2053

(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). 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 heart muscle. The number 1-4 distinguishes between the 4 different species.



Species 1-4

1 = Mouse

2 = Rat

3 = Rabbit

4 = Sheep

HM= Heart muscle

LU= Lung

LI= Liver

SP= Spleen

KI= Kidney

CN= Central nervous system

MS= Muscle, skeletal

BM= Bone marrow

TO= Tongue/Oral cavity

SK= Skin

ST= Stomach

SB= Small bowel

CO= Colon

OT= Ovary/Testis

UB= Urinary bladder

M*: Marking tissue (heart muscle tissue) for orientation of the MaxArray.

www.invitrogen.com

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

PI 75-2053

(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). 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.