

■ General Information

Applications

- Immunohistochemistry
 - TUNEL for apoptosis
- In situ hybridization (ISH)
 - mRNA
 - miRNA
 - Fluorescent In situ hybridization (FISH)

Storage and stability

- Individual slide is put in an air-tight pack with inert gas.
- If the slides are stored at 4C, they are good for up to one year.

How processed

- Tissues were initially fixed with formalin except for some of the animal tissues
- Then, dehydrated with gradient ethanol; typically 1 hour each progressive steps; 70%, 90%, 95%, 99%, 100% x 3 times.
- Cleared by xylene, three changes for 1 hour each.
- Infiltrated with 60°C paraffin, three changes for 1 hour each
- Sectioned by microtome in 4 µm thickness

Before use

- Dry slides for 1 hour in a oven at 60C.
- Dewax slides in xylene for 4 minutes x 5 times.
- Hydrate slides in 100%, 95% and 75% ethanol for 3 minutes x 2 times each.
- Immerse slides in tap water for 5 minutes.

Slide orientation

- In most of the slides with 59 or 60 cores, the orientation is as below unless indicated otherwise. #60 location is usually filled with carbon for orientation.

| | | | | | | | | | | |
|-------------|----|----|----|----|----|----|----|----|----|----|
| Shaded area | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 |
| | 11 | 12 | 13 | 14 | 15 | 16 | 17 | 18 | 19 | 20 |
| | 21 | 22 | 23 | 24 | 25 | 26 | 27 | 28 | 29 | 30 |
| | 30 | 32 | 33 | 34 | 35 | 36 | 37 | 38 | 39 | 40 |
| | 41 | 42 | 43 | 44 | 45 | 46 | 47 | 48 | 49 | 50 |
| | 51 | 52 | 53 | 54 | 55 | 56 | 57 | 58 | 59 | 60 |

■ Tissue types*

The "tissue type" column in the data sheet denotes the following categories.

1. normal tissue from a non-cancer patient
2. normal tissue from a cancer patient, but the cancer involves unrelated organ
3. normal tissue adjacent to the cancer
4. benign tumor
5. tumor of borderline malignancy or uncertain malignant potential
6. cancer

NBP2-30225 - Mouse Multi-Organ Tissue MicroArray (Normal)

| No. | Sex | Organ | Abnormal finding |
|-----|-----|----------------------------------|------------------|
| 1 | M | Skin | . |
| 2 | F | Skin | . |
| 3 | M | Spleen | . |
| 4 | F | Spleen | . |
| 5 | M | Skeletal muscle (abdominal wall) | . |
| 6 | F | Skeletal muscle (abdominal wall) | . |
| 7 | M | Trachea | . |
| 8 | F | Trachea | . |
| 9 | M | Lung | . |
| 10 | F | Lung | . |
| 11 | M | Heart | . |
| 12 | F | Heart | . |
| 13 | M | Salivary gland | . |
| 14 | F | Salivary gland | . |
| 15 | M | Liver | . |
| 16 | F | Liver | . |
| 17 | M | Pancreas | . |
| 18 | F | Pancreas | . |
| 19 | M | Esophagus | . |
| 20 | F | Esophagus | . |
| 21 | M | Forestomach | . |
| 22 | F | Forestomach | . |
| 23 | M | Glandular stomach | . |
| 24 | F | Glandular stomach | . |
| 25 | M | Duodenum | . |
| 26 | F | Duodenum | . |
| 27 | M | Jejunum | . |
| 28 | F | Jejunum | . |
| 29 | M | Ileum | . |
| 30 | F | Ileum | . |
| 31 | M | Colon | . |
| 32 | F | Colon | . |
| 33 | M | Cecum | . |
| 34 | F | Cecum | . |
| 35 | M | Rectum | . |
| 36 | F | Rectum | . |
| 37 | M | Kidney (cortex) | . |
| 38 | F | Kidney (cortex) | . |
| 39 | M | Kidney (medulla) | . |
| 40 | F | Kidney (medulla) | . |
| 41 | M | Bladder | . |
| 42 | F | Bladder | . |
| 43 | M | Prostate | . |
| 44 | M | Seminal vesicle | . |
| 45 | M | Testis | . |
| 46 | F | Uterus | . |
| 47 | F | Ovary | . |
| 48 | M | Adrenal gland | . |
| 49 | F | Adrenal gland | . |
| 50 | F | Thymus | . |
| 51 | M | Cerebrum | . |
| 52 | F | Cerebrum | . |
| 53 | M | Cerebellum | . |
| 54 | F | Cerebellum | . |
| 55 | M | Brain stem | . |
| 56 | M | Eyeball | . |
| 57 | F | Eyeball | . |
| 58 | M | Ear lobe | . |
| 59 | F | Ear lobe | . |
| 60 | . | Carbon | . |

ICR mice, 8 weeks old
 Perfusion fixation with 4% paraformaldehyde
 Suitable for in situ hybridization and immunohistochemistry