

■ 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

NBP2-30233 - Human Multi-tissue Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	Tissue type*
1	35	F	Skin	normal	3
2	46	F	Breast	normal	3
3	53	F	Spleen	normal	2
4	50	F	Lymph node	normal	2
5	57	M	Skeletal muscle	normal	2
6	39	M	Lung	normal	3
7	67	M	Heart	normal	2
8	69	M	Salivary gland, sublingual	normal	2
9	70	M	Liver	normal	1
10	65	M	Gallbladder	normal	3
11	73	F	Pancreas	normal	1
12	22	M	Tonsil	normal	3
13	76	F	Esophagus	normal	3
14	63	M	Stomach, antrum	normal	3
15	56	M	Stomach, fundus	normal	3
16	40	F	stomach, muscle layer	normal	3
17	44	M	Small bowel	normal	3
18	62	M	appendix	normal	3
19	73	M	Colon	normal	3
20	64	M	Rectum	normal	3
21	49	M	Kidney, cortex	normal	2
22	54	M	Kidney, medulla	normal	2
23	82	M	Urinary bladder	normal	1
24	32	M	Prostate	normal	1
25	70	M	Testis	normal	1
26	29	F	Endometrium	normal	1
27	47	F	Myometrium	normal	1
28	34	F	Placenta	normal	3
29	44	M	Adrenal gland	normal	1
30	39	M	Thyroid	normal	1
31	59	M	Cerebrum	normal	6
32	1	M	Cerebellum	normal	6
33	65	M	Skin	squamous cell carcinoma	6
34	36	M	Subcutis	liposarcoma	6
35	58	F	Breast	infiltrating duct carcinoma	6
36	34	F	Lymph node	Hodgkin lymphoma	6
37	54	F	Bone	osteosarcoma	6
38	61	M	Lung	adenocarcinoma	6
39	72	M	Lung	squamous cell carcinoma	6
40	41	F	Liver	cholangiocarcinoma	6
41	54	M	Liver	hepatocellular carcinoma	6
42	52	M	Liver	metastatic adenocarcinoma (from rectum)	6
43	77	M	Esophagus	squamous cell carcinoma	6
44	65	F	Stomach	adenocarcinoma	6
45	53	M	Stomach	malignant lymphoma, diffuse large B cell	6
46	40	F	Stomach	signet ring cell carcinoma	6
47	61	M	Duodenum	gastrointestinal stromal tumor, malignant	6
48	62	M	Descending colon	adenocarcinoma	6
49	73	M	Rectum	adenocarcinoma	6
50	57	M	Kidney	renal cell carcinoma	6
51	65	M	Urinary bladder	invasive urothelial carcinoma	6
52	63	M	Prostate	adenocarcinoma	6
53	35	M	Testis	seminoma	6
54	65	F	Uterine cervix	squamous cell carcinoma	6
55	69	F	Endometrium	adenocarcinoma	6
56	44	F	Ovary	metastatic adenocarcinoma (from stomach)	6
57	15	F	Ovary	mucinous cystadenocarcinoma	6
58	44	F	Ovary	serous cystadenoma of low malignant potential	6
59	69	F	Thyroid	papillary carcinoma	6
60	.	.	Carbon	.	.

uk: unknown

Tissue type*: see General information sheet

Supplementary data sheet contains survival data (available by web)