

## ■ 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
	31	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

## NBP2-30332 - Human Various Tissue MicroArray (Sarcoma)

No.	Sex	Age	Organ	Diagnosis	Tissue type*
1	F	50	Kidney	leiomyosarcoma	6
2	F	86	Soft tissue, knee	malignant fibrous histiocytoma	6
3	M	59	Soft tissue, chest wall	synovial sarcoma, biphasic type	6
4	F	58	Soft tissue, back	extraskelatal osteosarcoma	6
5	M	44	Retroperitoneum	liposarcoma, myxoid and round cell type	6
6	F	27	Uterus	leiomyosarcoma	6
7	M	63	Soft tissue, knee	malignant fibrous histiocytoma	6
8	M	66	Soft tissue and femur, thigh	malignant fibrous histiocytoma	6
9	F	27	Soft tissue, thigh	synovial sarcoma, monophasic type	6
10	M	50	Abdominal cavity	primitive neuroectodermal tumor	6
11	M	62	Soft tissue, thigh	malignant fibrous histiocytoma	6
12	M	62	Retroperitoneum	embryonal rhabdomyosarcoma	6
13	F	39	Abdominal cavity	leiomyosarcoma	6
14	M	19	Soft tissue, thigh	synovial sarcoma, biphasic type	6
15	M	46	Soft tissue, forearm	malignant peripheral nerve sheath tumor	6
16	M	41	Abdominal cavity	liposarcoma, myxoid type	6
17	M	32	Soft tissue, thigh	extraskelatal osteosarcoma	6
18	F	81	Pelvic cavity	leiomyosarcoma	6
19	F	54	Coccyx	chordoma	6
20	F	58	Retroperitoneum	leiomyosarcoma, epithelioid type	6
21	M	32	Soft tissue, hip	malignant mesenchymoma, metastatic	6
22	F	47	Retroperitoneum	liposarcoma, myxoid	6
23	M	38	Mesentery	liposarcoma, well differentiated	6
24	M	54	Soft tissue, thigh	malignant fibrous histiocytoma	6
25	F	54	Soft tissue, vertebra	malignant fibrous histiocytoma, metastatic	6
26	M	68	Soft tissue, thigh	leiomyosarcoma	6
27	M	15	Soft tissue, thigh	synovial sarcoma, biphasic type	6
28	F	13	Soft tissue, pelvis	primitive neuroectodermal tumor	6
29	M	31	Soft tissue, thigh	synovial sarcoma, biphasic type	6
30	F	22	Coccyx	chordoma	6
31	M	66	Soft tissue, arm	embryonal rhabdomyosarcoma	6
32	F	57	Soft tissue, face	dermatofibrosarcoma protuberans	6
33	F	67	Soft tissue, knee	malignant fibrous histiocytoma	6
34	M	67	Soft tissue, upper arm	malignant fibrous histiocytoma	6
35	F	53	Soft tissue, neck	malignant fibrous histiocytoma	6
36	M	48	Soft tissue, chest	malignant fibrous histiocytoma	6
37	F	30	Soft tissue, neck	malignant fibrous histiocytoma	6
38	M	36	Soft tissue, lower leg	round cell liposarcoma	6
39	F	79	Soft tissue, knee	malignant fibrous histiocytoma	6
40	M	42	Soft tissue, chest	leiomyosarcoma, epithelioid	6
41	M	43	Soft tissue, pelvis	synovial sarcoma, monophasic fibrous type	6
42	F	46	Soft tissue, thigh	leiomyosarcoma	6
43	F	67	Sacrum	chordoma	6
44	M	32	Pelvic cavity	primitive neuroectodermal tumor	6
45	F	65	Soft tissue, ankle	angiosarcoma	6
46	M	64	Soft tissue, thigh	liposarcoma	6
47	F	10	Soft tissue, cheek	alveolar rhabdomyosarcoma	6
48	F	71	Soft tissue, knee	pleomorphic rhabdomyosarcoma	6
49	M	46	Soft tissue, shoulder	synovial sarcoma	6
50	F	38	Small bowel	desmoplastic small round cell tumor	6
51	M	28	Soft tissue, thigh	epithelioid sarcoma	6
52	M	23	Soft tissue, thigh	fibrosarcoma	6
53	F	13	Pelvic cavity	neoplasm with perivascular epithelioid cell differentiation (PEComa)	6
54	M	32	Kidney	primitive neuroectodermal tumor	6
55	F	26	Femur	malignant giant cell tumor	6
56	M	56	Spermatic cord	leiomyosarcoma	6
57	M	44	Descending colon	malignant peripheral nerve sheath tumor	6
58	F	33	Abdominal wall	synovial sarcoma	6
59	F	49	Soft tissue, axilla	liposarcoma, well differentiated	6
60	.	.	Carbon	.	.