

■ 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

NBP2-30169 - Human Prostate Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	Gleason score	pTNM	Stage	Tissue type*
1	60	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
2	64	M	Prostate	adenocarcinoma	7	T2cNOMO	II B	6
3	71	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
4	64	M	Prostate	adenocarcinoma	10	T3aNOMO	III	6
5	59	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
6	65	M	Prostate	adenocarcinoma	8	T3bNOMO	III	6
7	73	M	Prostate	adenocarcinoma	7	T2cNOMO	II B	6
8	69	M	Prostate	adenocarcinoma	7	T2cNOMO	II B	6
9	62	M	Prostate	adenocarcinoma	7	T2cNOMO	II B	6
10	66	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
11	60	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
12	66	M	Prostate	adenocarcinoma	7	T3aNOMO	III	6
13	70	M	Prostate	adenocarcinoma	7	T3bNOMO	III	6
14	65	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
15	67	M	Prostate	adenocarcinoma	9	T3bN1MO	III	6
16	69	M	Prostate	adenocarcinoma	7	T3bNOMO	III	6
17	63	M	Prostate	adenocarcinoma	9	T3aN1MO	III	6
18	69	M	Prostate	adenocarcinoma	7	T3aNOMO	III	6
19	70	M	Prostate	adenocarcinoma	7	T3aNOMO	III	6
20	58	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
21	58	M	Prostate	adenocarcinoma	7	T3bNOMO	III	6
22	71	M	Prostate	adenocarcinoma	7	T2cNOMO	II B	6
23	70	M	Prostate	adenocarcinoma	7	T3bNOMO	III	6
24	59	M	Prostate	adenocarcinoma	6	T2bNOMO	II A	6
25	63	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
26	72	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
27	66	M	Prostate	adenocarcinoma	8	T3bNOMO	III	6
28	70	M	Prostate	adenocarcinoma	6	T3bNOMO	III	6
29	70	M	Prostate	adenocarcinoma	7	T2cNOMO	II B	6
30	68	M	Prostate	adenocarcinoma	8	T3bNOMO	III	6
31	63	M	Prostate	adenocarcinoma	10	T3bNOMO	III	6
32	57	M	Prostate	adenocarcinoma	7	T3bNOMO	III	6
33	72	M	Prostate	adenocarcinoma	8	T2cNOMO	II B	6
34	70	M	Prostate	adenocarcinoma	8	T3bNOMO	III	6
35	75	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
36	62	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
37	63	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
38	53	M	Prostate	adenocarcinoma	9	T3bNOMO	III	6
39	63	M	Prostate	adenocarcinoma	8	T3bNOMO	III	6
40	44	M	Prostate	adenocarcinoma	7	T3bNOMO	III	6
41	69	M	Prostate	normal (match of #8)	.	.	.	3
42	62	M	Prostate	normal (match of #9)	.	.	.	3
43	66	M	Prostate	normal (match of #12)	.	.	.	3
44	65	M	Prostate	normal (match of #14)	.	.	.	3
45	69	M	Prostate	normal (match of #18)	.	.	.	3
46	70	M	Prostate	normal (match of #19)	.	.	.	3
47	70	M	Prostate	normal (match of #29)	.	.	.	3
48	63	M	Prostate	normal (match of #31)	.	.	.	3
49	44	M	Prostate	normal (match of #40)	.	.	.	3
50	.	.	Carbon

TNM and Stage: AJCC Cancer Staging Manual (7th Edition)