

## ■ 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

### NBP3-11846 - Human Bladder Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	LN*	Tissue type **
1	UK	UK	Urinary bladder	Urothelial carcinoma	T4aN0M0	IIIA	0/0	6
2	43	M	Urinary bladder	Mucinous adenocarcinoma, urachal origin	T2bN0M0	II	0/0	6
3	80	M	Urinary bladder	Urothelial carcinoma	T2bN0M0	II	0/0	6
4	52	M	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/0	6
5	67	M	Urethra	Urothelial carcinoma	T2N0M0	II	0/0	6
6	71	M	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/0	6
7	74	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/0	6
8	52	M	Urinary bladder	Mucinous adenocarcinoma	T3aN0M0	IIIA	0/0	6
9	59	M	Urinary bladder	Urothelial carcinoma	T2bN0M0	II	0/0	6
10	66	M	Urinary bladder	Squamous cell carcinoma	T4aN1M0	IIIA	1/7	6
11	83	F	Urinary bladder	Urothelial carcinoma	T3aN0M0	IIIA	0/0	6
12	74	M	Urinary bladder	Urothelial carcinoma	T2bN0M0	II	0/12	6
13	68	F	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/1	6
14	63	M	Urinary bladder	Urothelial carcinoma	T3bN2M0	IIIB	3/19	6
15	64	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/0	6
16	34	M	Urinary bladder	Inflammatory myofibroblastic tumor	(4.5cm)	-	0/0	6
17	53	M	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/13	6
18	61	M	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/9	6
19	69	M	Urinary bladder	Urothelial carcinoma	T3bN2M0	IIIB	3/14	6
20	80	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/11	6
21	64	F	Urinary bladder	Urothelial carcinoma	T4aN0M0	IIIA	0/26	6
22	92	M	Urinary bladder	Urothelial carcinoma	T4aN0M0	IIIA	0/11	6
23	84	M	Urinary bladder	Urothelial carcinoma	T3aN0M0	IIIA	0/1	6
24	78	M	Urinary bladder	Urothelial carcinoma	T4aN0M0	IIIA	0/0	6
25	77	M	Urinary bladder	Urothelial carcinoma	T2bN0M0	II	0/13	6
26	71	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/14	6
27	64	M	Urinary bladder	Urothelial carcinoma	T2aN0M0	II	0/51	6
28	67	F	Urinary bladder	Urothelial carcinoma	T4aN0M0	IIIA	0/10	6
29	70	M	Urinary bladder	Urothelial carcinoma	T4aN0M0	IIIA	0/0	6
30	54	M	Urinary bladder	Urothelial carcinoma	T2bN0M0	II	0/0	6
31	77	M	Urinary bladder	Urothelial carcinoma	T4aN0M0	IIIA	0/0	6
32	65	M	Urinary bladder	Urothelial carcinoma	T4aN0M0	IIIA	0/37	6
33	70	M	Urinary bladder	Urothelial carcinoma	T4bN0M0	IVA	0/16	6
34	55	F	Urinary bladder	Urothelial carcinoma	T2aN0M0	II	0/34	6
35	58	M	Urinary bladder	Urothelial carcinoma	T2aN0M0	II	0/48	6
36	81	M	Urinary bladder	Urothelial carcinoma	T2aN0M0	II	0/24	6
37	72	F	Urinary bladder	Urothelial carcinoma	T4bN1M0	IVA	1/5	6
38	73	M	Urinary bladder	Urothelial carcinoma	T3bN2M0	IIIB	2/51	6
39	66	M	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/5	6
40	79	M	Urinary bladder	Urothelial carcinoma	T3bN2M0	IIIB	2/14	6
41	74	M	Urinary bladder	Urothelial carcinoma	T3bN1M0	IIIA	1/5	6
42	70	M	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/25	6
43	59	M	Urinary bladder	Urothelial carcinoma	TisN0M0	0is	0/27	6
44	49	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/10	6
45	93	M	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/0	6
46	69	F	Urinary bladder	Urothelial carcinoma	T3bN0M0	IIIA	0/40	6
47	59	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/18	6
48	55	M	Urinary bladder	Urothelial carcinoma	T3bN1M0	IIIA	1/28	6
49	58	F	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/15	6
50	65	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/37	6
51	78	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/0	6
52	52	F	Urinary bladder	Urothelial carcinoma	T2aN0M0	II	0/11	6
53	48	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/16	6
54	70	M	Urinary bladder	Squamous cell carcinoma	T3bN0M0	IIIA	0/17	6
55	73	M	Urinary bladder	Urothelial carcinoma	T3bN2M0	IIIB	12/35	6
56	61	M	Urinary bladder	Urothelial carcinoma	T4aN2M0	IIIB	11/33	6
57	70	M	Urinary bladder	Urothelial carcinoma	T1N0M0	I	0/23	6
58	65	M	Urinary bladder	Urothelial carcinoma	rT3aN0M0	IIIA	0/32	6
59	59	M	Urinary bladder	Urothelial carcinoma	T2bN0M0	II	0/25	6
60	.	.	Carbon	.	.	.	.	.

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