

■ 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

NBP2-30262 - Human Multi-tissue Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	NBP2-30215#	Tissue type*
1	56	M	Stomach	adenocarcinoma, poorly differentiated	T3N0M0	II	101	6
2	52	M	Stomach	adenocarcinoma, poorly differentiated	T3N1M0	IIIA	102	6
3	39	M	Stomach	adenocarcinoma, poorly differentiated	T3N2M0	IIIB	103	6
4	44	M	Stomach	adenocarcinoma, poorly differentiated	T3N3M0	IV	104	6
5	65	F	Stomach	mucinous adenocarcinoma	T3N1M0	IIIA	105	6
6	52	M	Stomach	adenocarcinoma, moderately differentiated	T3N1M0	IIIA	106	6
7	66	M	Stomach	adenocarcinoma, poorly differentiated	T3N2M0	IIIB	107	6
8	61	M	Stomach	adenocarcinoma, poorly differentiated	T3N1M0	IIIA	108	6
9	39	F	Stomach	signet ring cell carcinoma	T3N0M0	II	109	6
10	60	F	Stomach	adenocarcinoma, moderately differentiated	T3N1M0	IIIA	110	6
11	64	M	Esophagus	adenocarcinoma, poorly differentiated	T3N0M0	IIA	111	6
12	60	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III	112	6
13	73	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III	113	6
14	52	F	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III	114	6
15	77	M	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	IIA	115	6
16	63	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III	116	6
17	47	M	Esophagus	squamous cell carcinoma, well differentiated	T2N1M0	IIB	117	6
18	58	M	Esophagus	basaloid squamous cell carcinoma	T3N0M0	IIA	118	6
19	60	F	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	IIA	119	6
20	69	M	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	IIA	120	6
21	69	M	Lung	adenocarcinoma, well differentiated	T2N0M0	IB	121	6
22	71	M	Lung	large cell carcinoma	T2N0M0	IB	122	6
23	65	M	Lung	squamous cell carcinoma, moderately differentiated	T2N0M0	IB	123	6
24	62	M	Lung	squamous cell carcinoma, moderately differentiated	T2N1M0	IIB	124	6
25	69	M	Lung	squamous cell carcinoma, moderately differentiated	T4N0M0	IIIB	125	6
26	58	M	Lung	adenocarcinoma, well differentiated	T2N0M0	IB	126	6
27	58	M	Lung	squamous cell carcinoma, moderately differentiated	T2N1M0	IIB	127	6
28	61	M	Lung	squamous cell carcinoma, moderately differentiated	T3N2M0	IIIA	128	6
29	52	M	Lung	squamous cell carcinoma, well differentiated	T2N1M0	IIB	129	6
30	56	M	Lung	squamous cell carcinoma, well differentiated	T2N0M0	IB	130	6
31	66	M	Sigmoid colon	adenocarcinoma, well differentiated	T3N0M0	IIA	131	6
32	50	M	Rectum	adenocarcinoma, moderately differentiated	T3N0M0	IIA	132	6
33	40	M	Cecum	adenocarcinoma, moderately differentiated	T3N0M0	IIA	133	6
34	45	M	Ascending colon	adenocarcinoma, well differentiated	T3N0M0	IIA	134	6
35	43	F	Rectum	mucinous adenocarcinoma	T3N2M0	IIIC	135	6
36	44	F	Ascending colon	adenocarcinoma, moderately differentiated	T3N1M0	IIIA	136	6
37	52	F	Rectum	adenocarcinoma, moderately differentiated	T3N0M0	IIA	137	6
38	49	M	Ascending colon	adenocarcinoma, moderately differentiated	T3N1M0	IIIA	138	6
39	61	F	Rectum	adenocarcinoma, moderately differentiated	T3N1M0	IIIA	139	6
40	52	F	Rectum	adenocarcinoma, moderately differentiated	T2N0M0	I	140	6
41	29	F	Thyroid	papillary carcinoma	T2N0M0	I	.	6
42	20	F	Thyroid	papillary carcinoma	T2N0M0	I	142	6
43	15	F	Thyroid	papillary carcinoma	T2N0M0	I	143	6
44	60	F	Thyroid	papillary carcinoma	T2N0M0	II	144	6
45	47	F	Thyroid	papillary carcinoma	T2N0M0	II	145	6
46	48	F	Thyroid	papillary carcinoma	T3N1aM0	III	146	6
47	43	F	Thyroid	papillary carcinoma	T1N0M0	I	147	6
48	56	F	Thyroid	papillary carcinoma	T2N1bM0	IVA	148	6
49	55	F	Thyroid	papillary carcinoma	T3N0M0	III	149	6
50	34	F	Thyroid	papillary carcinoma	T3N0M0	I	150	6
51	57	F	Kidney	renal cell carcinoma, clear cell type	T1aN0M0	I	151	6
52	50	F	Kidney	renal cell carcinoma, clear cell type	T2N0M0	II	152	6
53	27	M	Kidney	nephrogenic carcinoma	T2N0M0	II	153	6
54	56	M	Kidney	renal cell carcinoma, clear cell type	T1aN0M0	I	154	6
55	47	M	Kidney	renal cell carcinoma, granular cell type	T2N0M0	II	155	6
56	63	M	Kidney	renal cell carcinoma	T2N0M0	II	156	6
57	53	M	Kidney	renal cell carcinoma, clear cell type	T3aN0M0	III	157	6
58	62	M	Kidney	renal cell carcinoma, clear cell type	T2N0M0	II	158	6
59	67	F	Kidney	renal cell carcinoma, clear cell type	T3aN0M0	III	.	6
60	.	.	Carbon

#: The normal tissue in NBP2-30215 of corresponding number is from the identical patient. TNM and Stage: AJCC Cancer Staging Manual (6th Edition)