

■ General Information of the Tissue Array Slides

Applications

- Immunohistochemistry
- Fluorescent in situ hybridization (FISH)
- mRNA in situ hybridization
- miRNA in situ hybridization
- TUNEL for apoptosis
- Nucleic acid extraction

Storage and stability

- Individual slide is put in an air-tight pack with inert gas.
- Tissue array slides are shipped at room temperature.
- If the slides are stored at 4 degree, 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 and put on New Silane III slide by Muto Pure Chemicals.

Before use

- Dry slides for 1 hour in a oven at 60 degree.
- 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

NBP2-30308 - Human Stomach Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	No. of CQN3*	Tissue type**
1	57	M	Stomach	adenocarcinoma, poorly differentiated	T4bN2M0	IIIC	101	6
2	59	M	Stomach	adenocarcinoma, poorly differentiated	T3N2M0	IIIA	102	6
3	64	M	Stomach	adenocarcinoma, moderately differentiated	T1bN2M0	IIA	103	6
4	45	M	Stomach	adenocarcinoma, poorly differentiated	T3N3bM0	IIIB	104	6
5	67	M	Stomach	adenocarcinoma, poorly differentiated	T3N1M0	IIB	105	6
6	50	M	Stomach	adenocarcinoma, moderately differentiated	T3N3aM0	IIIB	106	6
7	67	M	Stomach	adenocarcinoma, moderately differentiated	T4aN3bM0	IIIC	107	6
8	55	F	Stomach	adenocarcinoma, poorly differentiated	T2N0M0	IB	108	6
9	51	F	Stomach	signet ring cell carcinoma	T4aN3aM0	IIIC	109	6
10	65	M	Stomach	adenocarcinoma, poorly differentiated	T4aN0M0	IIB	110	6
11	70	M	Stomach	adenocarcinoma, moderately differentiated	T4aN2M0	IIIB	111	6
12	46	M	Stomach	adenocarcinoma, poorly differentiated	T3N3aM0	IIIB	112	6
13	49	M	Stomach	adenocarcinoma, moderately differentiated	T3N2M0	IIIA	113	6
14	52	M	Stomach	papillary adenocarcinoma	T2N0M0	IB	114	6
15	62	M	Stomach	signet ring cell carcinoma	T3N3aM0	IIIB	115	6
16	66	M	Stomach	adenocarcinoma, poorly differentiated	T4aN3bM0	IIIC	116	6
17	55	M	Stomach	adenocarcinoma, moderately differentiated	T3N3aM1	IV	117	6
18	51	F	Stomach	signet ring cell carcinoma	T3N3aM0	IIIB	118	6
19	60	M	Stomach	adenocarcinoma, moderately differentiated	T3N3aM0	IIIB	119	6
20	64	M	Stomach	adenocarcinoma, moderately differentiated	T4aN3bM1	IV	120	6
21	53	M	Stomach	signet ring cell carcinoma	T2N0M0	IB	121	6
22	64	M	Stomach	mucinous adenocarcinoma	T3N1M0	IIB	122	6
23	71	M	Stomach	undifferentiated carcinoma	T3N0M0	IIA	123	6
24	69	M	Stomach	adenocarcinoma, poorly differentiated	T4aN3aM0	IIIC	124	6
25	64	M	Stomach	adenocarcinoma, moderately differentiated	T4aN3aM0	IIIC	125	6
26	63	M	Stomach	signet ring cell carcinoma	T4bN3aM0	IIIC	126	6
27	39	F	Stomach	adenocarcinoma, moderately differentiated	T3N3bM0	IIIB	127	6
28	71	M	Stomach	adenocarcinoma, moderately differentiated	T3N3bM0	IIIB	128	6
29	63	M	Stomach	adenocarcinoma, moderately differentiated	T4aN3aM0	IIIC	129	6
30	77	M	Stomach	adenocarcinoma, moderately differentiated	T3N3bM0	IIIB	130	6
31	44	F	Stomach	adenocarcinoma, poorly differentiated	T4aN3bM1	IV	131	6
32	59	M	Stomach	adenocarcinoma, poorly differentiated	T3aN0M0	IIA	132	6
33	42	M	Stomach	adenocarcinoma, moderately differentiated	T3N3aM0	IIIB	133	6
34	31	F	Stomach	adenocarcinoma, poorly differentiated	T4aN1M1	IV	134	6
35	54	F	Stomach	adenocarcinoma, poorly differentiated	T2N2M0	IIB	135	6
36	65	M	Stomach	adenocarcinoma, poorly differentiated	T2N1M0	IIA	136	6
37	34	F	Stomach	adenocarcinoma, poorly differentiated	T3N0M0	IIA	137	6
38	74	M	Stomach	adenocarcinoma, poorly differentiated	T4aN3bM0	IIIC	138	6
39	63	M	Stomach	adenocarcinoma, moderately differentiated	T4aN3bM0	IIIC	139	6
40	49	M	Stomach	adenocarcinoma, poorly differentiated	T4aN3aM0	IIIC	140	6
41	48	M	Stomach	adenocarcinoma, poorly differentiated	T4aN3bM0	IIIC	141	6
42	52	F	Stomach	adenocarcinoma, moderately differentiated	T3N2M0	IIIA	142	6
43	49	F	Stomach	signet ring cell carcinoma	T4aN2M0	IIIB	143	6
44	76	M	Stomach	adenocarcinoma, poorly differentiated	T2N0M0	IB	144	6
45	55	M	Stomach	adenocarcinoma, well differentiated	T1aN0M0	IA	145	6
46	43	M	Stomach	adenocarcinoma, moderately differentiated	T4aN2M0	IIIB	146	6
47	59	M	Stomach	adenocarcinoma, moderately differentiated	T4aN2M0	IIIB	147	6
48	75	M	Stomach	adenocarcinoma, moderately differentiated	T4aN1M0	IIIA	148	6
49	62	F	Stomach	adenocarcinoma, poorly differentiated	T1bN0M0	IA	149	6
50	72	F	Stomach	adenocarcinoma, poorly differentiated	T4aN0M0	IIB	150	6
51	56	M	Stomach	adenocarcinoma, poorly differentiated	T2N1M0	IIA	151	6
52	47	M	Stomach	adenocarcinoma, moderately differentiated	T3N1M0	IIB	152	6
53	59	F	Stomach	adenocarcinoma, well differentiated	T4aN3bM1	IV	153	6
54	42	M	Stomach	adenocarcinoma, poorly differentiated	T3N2M0	IIIA	154	6
55	46	M	Stomach	adenocarcinoma, moderately differentiated	T3N1M0	IIB	155	6
56	68	M	Stomach	adenocarcinoma, well differentiated	T1bN1M0	IB	156	6
57	64	F	Stomach	adenocarcinoma, poorly differentiated	T4aN3aM0	IIIC	157	6
58	64	F	Stomach	adenocarcinoma, moderately differentiated	T3N3bM0	IIIB	158	6
59	60	M	Stomach	adenocarcinoma, well differentiated	T2N0M0	IB	159	6
60	.	.	Carbon

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