

■ 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

NBP2-30219 - Human Gastrointestinal Tissue MicroArray (Stromal Tumor)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	Months*	Tissue type**
1	45	F	Jejunum	GIST, high risk	T3N0M0	III B	.	6
2	57	M	Jejunum	GIST, high risk	T4N0M0	III B	.	6
3	77	F	Stomach	GIST, high risk	T4N0M0	III B	.	6
4	60	F	Stomach	GIST, high risk	T3N0M0	III A	.	6
5	59	M	Jejunum	GIST, high risk	T3N0M0	III B	.	6
6	58	M	Small intestine	GIST, high risk	T4N1M0	IV	.	6
7	65	F	Stomach	GIST, low risk	T3N0M0	I B	.	4
8	61	M	Stomach	GIST, high risk	T4N1M0	IV	.	6
9	65	F	Stomach	GIST, high risk	T4N0M0	II	.	6
10	44	M	Small intestine	GIST, high risk	T3N0M0	III B	.	6
11	53	F	Duodenum	GIST, intermediate risk	T3N0M0	II	.	5
12	50	M	Stomach	GIST, low risk	T2N0M0	I A	.	4
13	68	M	Stomach	GIST, high risk	T4N0M0	III B	.	6
14	34	M	Duodenum	GIST, high risk	T2N0M0	II	.	6
15	61	M	Stomach	GIST, intermediate risk	T3N0M0	I B	.	5
16	58	F	Stomach	GIST, high risk	T3N0M0	III A	.	6
17	57	F	Stomach	GIST, low risk	T2N0M0	I A	.	4
18	40	M	Stomach	GIST, high risk	T4N0M0	III B	.	6
19	66	F	Rectum	GIST, high risk	T3N0M0	III B	.	6
20	45	F	Abdominal cavity	GIST, high risk	T4N0M1	IV	.	6
21	66	M	Stomach	GIST, high risk	T4N0M0	III B	.	6
22	83	F	Stomach	GIST, high risk	T4N0M0	III B	.	6
23	60	F	Jejunum	GIST, intermediate risk	T3N0M0	II	.	5
24	65	M	Jejunum	GIST, high risk	T4N0M0	III B	.	6
25	56	F	Duodenum	GIST, low risk	T2N0M0	I	.	4
26	85	F	Stomach	GIST, high risk	T4N0M0	II	.	6
27	67	F	Duodenum	GIST, intermediate risk	T3N0M0	II	.	5
28	52	M	Stomach	GIST, high risk	T4N0M0	III B	.	6
29	65	M	Stomach	GIST, low risk	T2N0M0	I A	.	4
30	65	M	Ileum	GIST, intermediate risk	T3N0M0	II	.	5
31	45	F	Stomach	GIST, low risk	T2N0M0	I A	.	4
32	55	M	Rectum	GIST, high risk	T3N0M0	III B	.	6
33	71	M	Stomach	GIST, high risk	T4N0M0	II	.	6
34	67	F	Stomach	GIST, intermediate risk	T2N0M0	II	.	5
35	62	F	Stomach	GIST, intermediate risk	T3N0M0	I B	.	5
36	48	M	Stomach	GIST, high risk	T4N0M0	III B	.	6
37	80	M	Small intestine	GIST, low risk	T2N0M0	I	.	4
38	60	M	Small intestine	GIST, intermediate risk	T3N0M0	II	.	5
39	43	M	Duodenum	GIST, low risk	T3N0M0	II	.	4
40	53	F	Stomach	GIST, high risk	T3N0M0	III A	.	6
41	78	M	Liver	spindle cell sarcoma, metastatic	M1	IV	0	6
42	81	F	Lung	leiomyosarcoma, metastatic	M1	IV	0	6
43	65	M	Omentum	GIST, recurred	T4N0M0	III B	75	6
44	80	M	Liver	GIST, metastatic	M1	IV	48	6
45	64	M	Mesentery	GIST, metastatic	T4N0M0	III B	28	6
46	67	M	Mesentery	GIST, metastatic	T2N0M0	I	48	6
47	35	M	Ileum	GIST, recurred	T3N0M0	III B	36	6
48	62	M	Omentum	GIST, metastatic	T4N0M0	III B	0	6
49	45	F	Liver	GIST, metastatic (match of #1)	M1	IV	80	6
50	57	M	Mesentery	GIST, low risk, recurred (match of #2)	T2N0M0	I	16	4
51	77	F	Stomach	normal (match of #3)	.	.	.	3
52	60	F	Stomach	normal (match of #4)	.	.	.	3
53	59	M	Jejunum	normal (match of #5)	.	.	.	3
54	58	M	Small intestine	normal (match of #6)	.	.	.	3
55	61	M	Stomach	normal (match of #8)	.	.	.	3
56	44	M	Small intestine	normal (match of #10)	.	.	.	3
57	68	M	Stomach	normal (match of #13)	.	.	.	3
58	66	F	Rectum	normal (match of #19)	.	.	.	3
59	65	M	Jejunum	normal (match of #24)	.	.	.	3
60	.	.	Carbon

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

Months*: interval between primary and metastatic/recurrent cancer