

■ 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

NBP2-30256 - Human Colon Tissue MicroArray (Rectal Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	stage	No. of NBP2-30213#	Tissue type*
1	60	F	Ascending colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	101	6
2	77	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T2N0M0	I	102	6
3	51	M	Transverse colon	adenocarcinoma, poorly differentiated	T3N0M0	IIA	103	6
4	38	F	Transverse colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	104	6
5	56	M	Sigmoid colon	adenocarcinoma, well differentiated	T3N2aM0	IIIB	105	6
6	46	F	Cecum	adenocarcinoma, moderately differentiated	T2N1aM0	IIIA	106	6
7	61	M	Cecum	mucinous carcinoma	T3N0M0	IIA	107	6
8	70	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	108	6
9	36	F	Descending colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	109	6
10	46	F	Rectum	adenocarcinoma, well differentiated	T3N0M0	IIA	110	6
11	62	F	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	111	6
12	65	M	Ascending colon	adenocarcinoma, well differentiated	T3N2aM0	IIIB	112	6
13	65	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	113	6
14	72	F	Sigmoid colon	adenocarcinoma, well differentiated	T3N1bM0	IIIB	114	6
15	63	M	Ascending colon	adenocarcinoma, well differentiated	T3N0M0	IIA	115	6
16	64	M	Sigmoid colon	adenocarcinoma, well differentiated	T3N1aM0	IIIB	116	6
17	64	F	Rectum	adenocarcinoma, moderately differentiated	T2N0M0	I	117	6
18	71	F	Descending colon	mucinous adenocarcinoma	T3N0M0	IIA	118	6
19	57	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N0M1a	IVA	119	6
20	58	M	Cecum	adenocarcinoma, moderately differentiated	T3N0M0	IIA	120	6
21	56	M	Ascending colon	mucinous carcinoma	T3N0M0	IIA	121	6
22	57	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T4bN0M0	IIC	122	6
23	56	M	Ascending colon	adenocarcinoma, moderately differentiated	T3N1bM0	IIIB	123	6
24	36	M	Transverse colon	adenocarcinoma, moderately differentiated	T3N2bM0	IIIC	124	6
25	52	M	Rectum	adenocarcinoma, well differentiated	T4aN0M0	IIB	125	6
26	62	M	Ascending colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	126	6
27	57	F	Sigmoid colon	adenocarcinoma, well differentiated	T3N1bM0	IIIB	127	6
28	78	M	Sigmoid colon	adenocarcinoma, well differentiated	T3N1aM0	IIIB	128	6
29	71	M	Ascending colon	mucinous adenocarcinoma	T3N0M0	IIA	129	6
30	65	M	Ascending colon	adenocarcinoma, well differentiated	T3N1aM0	IIIB	130	6
31	61	M	Transverse colon	mucinous carcinoma	T3N0M0	IIA	131	6
32	52	M	Rectum	adenocarcinoma, moderately differentiated	T3N0M1a	IVA	132	6
33	52	F	Sigmoid colon	adenocarcinoma, moderately differentiated	T2N2aM0	IIIB	133	6
34	62	M	Descending colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	134	6
35	72	M	Transverse colon	adenocarcinoma, moderately differentiated	T3N2bM0	IIIC	135	6
36	54	M	Ascending colon	adenocarcinoma, poorly differentiated	T3N2bM0	IIIC	136	6
37	78	M	Ascending colon	adenocarcinoma, moderately differentiated	T3N1aM0	IIIB	137	6
38	64	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N1bM0	IIIB	138	6
39	59	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T4bN0M0	IIC	139	6
40	57	M	Ascending colon	adenocarcinoma, moderately differentiated	T3N0M1a	IVA	140	6
41	46	F	Ascending colon	adenocarcinoma, poorly differentiated	T3N2bM0	IIIC	141	6
42	64	M	Sigmoid colon	adenocarcinoma, well differentiated	T3N0M0	IIA	142	6
43	65	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T4bN2bM0	IIIC	143	6
44	76	F	Ascending colon	small cell carcinoma	T3N0M0	IIA	144	6
45	68	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N1aM0	IIIB	145	6
46	76	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N1bM1a	IVA	146	6
47	60	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	147	6
48	60	F	Rectum	adenocarcinoma, moderately differentiated	T4bN0M0	IIC	148	6
49	67	M	Descending colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	149	6
50	58	M	Ascending colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	150	6
51	57	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	151	6
52	55	M	Transverse colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	152	6
53	86	F	Ascending colon	adenocarcinoma, moderately differentiated	T3N2aM0	IIIB	153	6
54	58	M	Rectum	adenocarcinoma, moderately differentiated	T3N0M0	IIA	154	6
55	66	M	Ascending colon	mucinous adenocarcinoma	T3N0M0	IIA	155	6
56	35	M	Rectum	adenocarcinoma, moderately differentiated	T2N1aM0	IIIA	156	6
57	66	M	Sigmoid colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	157	6
58	60	M	Transverse colon	adenocarcinoma, well differentiated	T3N0M0	IIA	158	6
59	70	M	Descending colon	adenocarcinoma, moderately differentiated	T3N0M0	IIA	159	6
60	.	.	Carbon

#: The normal tissue in NBP2-30213 array of corresponding number is from the identical patient TNM and stage: AJCC Cancer Staging Manual (7th Edition)