

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 μ m 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.

	1	2	3	4	5	6	7	8	9	10
	11	12	13	14	15	16	17	18	19	20
Chadad area	21	22	23	24	25	26	27	28	29	30
Shaueu area	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

NBP2-30329 - Human Various Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	Tissue type
1	44	Μ	Oral cavity, tongue	squamous cell carcinoma, well differentiated	T2N0M0		6
2	55	M	Oral cavity, tongue	squamous cell carcinoma, well differentiated	T2N0M0		6
3	55 40	IVI M	Oral cavity, lower lip	squamous cell carcinoma, well differentiated	T2N2bM0	111	6
4	38	M	Oral cavity, marilla	squamous cell carcinoma, weir direrentiated	T3N1M0		6
6	45	M	Oral cavity, maxilla	squamous cell carcinoma, moderately differentiated	T2N0M0	11	6
7	58	M	Oral cavity, tonsil	squamous cell carcinoma, well differentiated	T3N0M0		6
8	23	Μ	Nasopharynx	angiofibroma	T1N0M0		4
9	26	М	Nasopharynx	olfactory neuroblastoma	T1N0M0	1	6
10	19	F	Nasopharynx	olfactory neuroblastoma	T1N0M0		6
11	35	M	Nasopharynx	olfactory neuroblastoma, recurrent, extension to orbit	T1N0M0		6
12	32 52	F	Nasopharyny	undifferentiated carcinoma	T1N0M0		6
14	62	M	Salivary gland parotid		T3N0M0		6
15	53	M	Salivary gland, parotid	acinic cell carcinoma	T3N1M0		6
16	56	F	Salivary gland, parotid	acinic cell carcinoma	T2N0M0		6
17	51	F	Salivary gland, parotid	adenoid cystic carcinoma	T4aN0M0	IVA	6
18	72	F	Salivary gland, submandibular	adenoid cystic carcinoma	T4aN0M0	IVA	6
19	53	F	Salivary gland	adenoid cystic carcinoma, metastasis to breast	M1	IVC	6
20	21	F	Salivary gland, parotid	mucoepidermoid carcinoma, low grade	T3N2bM0	IVA	6
21	51 61	Г	Salivary gland, parotid	enithelial myoenithelial carcinoma	T3N0M0		6
23	42	M	Salivary gland, parotid	myoepithelioma, malignant, recurrent	T2N0M0	11	6
24	32	F	Salivary gland, parotid	carcinoma ex pleomorphic adenoma	T2N0M0	11	6
25	35	F	Salivary gland, submandibular	carcinoma ex pleomorphic adenoma	T4aN0M0	IVA	6
26	64	М	Salivary gland, parotid	carcinoma ex pleomorphic adenoma	T1N0M0		6
27	80	F	Salivary gland, parotid	carcinoma ex pleomorphic adenoma	T3N0M0		6
28	57	M	Salivary gland, parotid	squamous cell carcinoma	T3N2bM0	IVA	6
29	52	IVI M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0		6
30	62	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0		6
32	54	F	Stomach	adenocarcinoma, well differentiated	T3N3M0	IV	6
33	52	M	Stomach	adenocarcinoma, well differentiated	T3N0M0	11	6
34	65	Μ	Stomach	adenocarcinoma, moderately differentiated	T3N0M0		6
35	50	Μ	Stomach	adenocarcinoma, moderately differentiated	T3N3M0	IV	6
36	61	Μ	Stomach	adenocarcinoma, moderately differentiated	T3N1M0	IIIA	6
37	60	M	Stomach	adenocarcinoma, moderately differentiated	T2aN0M0	I B	6
38	62	IVI M	Stomach	adenocarcinoma, moderately differentiated	T2N1M0		6
40	53	M	Stomach	adenocarcinoma, poorly differentiated	T3N3M0	IV	6
41	62	M	Stomach	adenocarcinoma, poorly differentiated	T2aN0M0	I B	6
42	55	F	Stomach	adenocarcinoma, poorly differentiated	T3N0M0		6
43	32	F	Stomach	adenocarcinoma, poorly differentiated	T3N1M0	IIIA	6
44	56	F	Stomach	adenocarcinoma, poorly differentiated	T3N2M0	IIIB	6
45			Carbon	· · · · · · · · · · · · · · · · · · ·			
46	52	IVI M	Stomach	adenocarcinoma, poorly differentiated	T2N2M0	IV	6
47	68	F	Stomach	adenocarcinoma, poorly differentiated	T2hN1M1	IV	6
49	55	M	Stomach	adenocarcinoma, poorly differentiated	T3N1M0	IIIA	6
50	52	M	Stomach	adenocarcinoma, poorly differentiated	T3N2M0	IIIB	6
51	58	Μ	Stomach	adenocarcinoma, mucinous	T3N1M0	IIIA	6
52	71	М	Stomach	adenocarcinoma, mucinous	T3N3M0	IV	6
53	43	F	Stomach	adenocarcinoma, metastasis to ovary	M1	IV	6
54	53	F	Stomach	adenocarcinoma, metastasis to ovary	M1	IV	6
55	4/ 29	F	Stomach	adenocarcinoma, metastasis to ovary	M1	IV IV	6
57	61	F	Stomach	adenocarcinoma, metastasis to ovary	M1	IV	6
58	36	F	Stomach	adenocarcinoma, metastasis to ovary, signet ring cell	M1	IV	6
59	39	F	Stomach	adenocarcinoma, metastasis to ovary, mucinous	M1	IV	6
60	37	F	Stomach	adenocarcinoma, metastasis to ovary, signet ring cell	M1	IV	6
61	50	Μ	Stomach	adenocarcinoma, metastasis to lymph node	T3N3M0	IV	6
62	41	F	Stomach	adenocarcinoma, metastasis to lymph nodes, papillary	T1N3M0	IV	6
63	55	F	Stomach	gastrointestinal stromal tumor	T1aN0M0		3
64	39	IVI M	Small Intestine, jejunum	gastrointestinal stromal tumor, malignant		0	6
66	55	M	Small intestine, jejunum	gasironitestinal stromal tumor, malignant	T2bN1M0	IV	6
67	42	M	Small intestine	gastrointestinal stromal tumor, malignant	T2bN0M0	0	6
68	61	M	Small intestine, duodenum	adenocarcinoma, metastasis to lymph node	T4N1M1	IV	6
69	44	F	Colon, sigmoid	villotubular adenomas	T1N0M0		4
70	52	М	Colon, ascending	adenocarcinoma, well differentiated	T3N0M0	II A	6
71	71	F	Colon, sigmoid	adenocarcinoma, well differentiated	T3N0M0	II A	6
72	82	F	Colon, cecum	adenocarcinoma, well differentiated	T3N2M0	IIIC	6
73	50	F	Colon, sigmoid	adenocarcinoma, Well differentiated			6
75	40 5/	F	Colon	adenocarcinoma, well differentiated	T3N1WU T3N1M1	III D IV	6
76	73	F	Colon	adenocarcinoma, well differentiated. recurrent			6
_77	62	F	Colon, descending	adenocarcinoma, moderately differentiated	T3N0M0	A	6
78	46	Μ	Colon, sigmoid	adenocarcinoma, moderately differentiated	T3N0M0	A	6
79	73	F	Colon, cecum	adenocarcinoma, moderately differentiated	T3N0M0	A	6
80	72	F	Colon, cecum	adenocarcinoma, moderately differentiated	T3N1M0	IIIB	6

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	Tissue type
81	37	F	Colon, ascending	adenocarcinoma, poorly differentiated	T3N1M0	III B	6
82	44	М	Colon, ascending	adenocarcinoma, mucinous	T3N2M0	III C	6
83	69	F	Colon, transverse	adenocarcinoma, mucinous	T3N0M0	IIA	6
85	42	F	Colon	adenocarcinoma, metastasis to uterus, mucinous	M1	IV	6
86	59	M		adenocarcinoma, metastasis to deeras, indenodas	M1	IV	6
87	51	F	Rectum	adenocarcinoma, well differentiated	T3N2M0	IIIC	6
88	75	F	Rectum	adenocarcinoma, well differentiated	T2N0M0		6
89	43	М	Rectum	adenocarcinoma, well differentiated	T3N1M0	IIIB	6
90	60	М	Rectum	adenocarcinoma, well differentiated	T3N1M0	IIIB	6
91	55	F	Rectum	adenocarcinoma, well differentiated	T3N1M0	IIIB	6
92	63	F	Rectum	adenocarcinoma, moderately differentiated	T 3N2IVIO		6
93	63	F	Rectum	adenocarcinoma, moderately differentiated	T3N0N0		6
95			Carbon		131101110		
96	70	M	Rectum	adenocarcinoma, moderately differentiated	T3N0M0	A	6
97	61	F	Anus	squamous cell carcinoma	T4N0M0	IIIA	6
98	63	М	Liver	cholangiocarcinoma, poorly differentiated	T1N0M0	Ι	6
99	41	М	Liver	cholangiocarcinoma, moderately differentiated	T4N0M0	III B	6
100	56	М	Liver	cholangiocarcinoma, moderately differentiated	T4N0M0	IIIB	6
101	56	M	Liver	hepatocellular carcinoma, well differentiated	T1N0M0	1	6
102	54	M	Liver	hepatocellular carcinoma, well differentiated	T2N0M0	11	6
103	62	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0		6
104	45	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	11	6
106	51	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0		6
107	77	М	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	П	6
108	59	F	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0		6
109	64	М	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0		6
110	67	F	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	Ш	6
111	53	М	Liver	hepatocellular carcinoma, metastasis to abdominal cavity	M1	IV	6
112	52	M	Liver	hepatocellular carcinoma, metastasis to abdominal wall	M1	IV	6
113	51	F	Liver Biliary, common bile duct	nepatocellular carcinoma, metastasis to pelvic cavity	MT T2N1M0	IV	6
114	67	F	Biliary, common blie duct	adenocarcinoma, well differentiated	T3N0M0		6
116	45	F	Biliary common bile duct	adenocarcinoma, moderately differentiated	T3N0M0		6
117	64	F	Biliary, gallbladder	adenocarcinoma, moderately differentiated	T3N1M0	B	6
118	52	F	Biliary, gallbladder	adenocarcinoma, poorly differentiated	T3N0M0	A	6
119	66	F	Biliary, gallbladder	adenosquamous cell carcinoma	T3N0M0	A	6
120	71	М	Pancreas	adenocarcinoma, moderately differentiated	T3N1M0	II B	6
121	54	М	Pancreas	adenocarcinoma, moderately differentiated	T3N1M0	II B	6
122	57	F	Pancreas	adenocarcinoma, moderately differentiated	T4N0M0		6
123	/1	E E	Pancreas	adenocarcinoma, moderately differentiated	T4N0M0	111	6
124	42	M		squamous cell carcinoma, well differentiated	T4N1W0	IV A	6
125	81	F	Larvnx	squamous cell carcinoma, well differentiated	T3N0M0		6
127	71	M	Larynx	squamous cell carcinoma, well differentiated	T4aN1M0	IVA	6
128	67	М	Larynx	squamous cell carcinoma, well differentiated	T4aN0M0	IVA	6
129	52	М	Larynx	squamous cell carcinoma, well differentiated	T1N2cM0	IVA	6
130	54	М	Larynx	squamous cell carcinoma, well differentiated	T3N0M0		6
131	60	M	Larynx	squamous cell carcinoma, well differentiated	T4aN0M0	IVA	6
132	66 57	IVI M	Larynx	squamous cell carcinoma, well differentiated	T J NUMU	111	6
133	46	M	Larynx	squamous cell carcinoma, well differentiated	T4aN2aM0	IV A	6
135	66	M		squamous cell carcinoma, well differentiated	T2N1M0		6
136	57	M	Larynx	squamous cell carcinoma, moderately differentiated	T4aN0M0	IVA	6
137	64	М	Larynx	squamous cell carcinoma, moderately differentiated	T4aN0M0	IVA	6
138	68	М	Larynx	squamous cell carcinoma, moderately differentiated	T4aN2bM0	IVA	6
139	57	М	Larynx	squamous cell carcinoma, metastasis to lung	M1	IVC	6
140	62	М	Lung	adenocarcinoma, well differentiated	T2N0M0	I B	6
141	26	M	Lung	adenocarcinoma, well differentiated	T2N1M0	IIB	6
142	5/	M	Lung	adenocarcinoma, well differentiated	T2N2M0	III A	6
143	0U 71	Г M		adenocarcinoma, modetately differentiated		III A	0
145		171	Carbon		121101010		0
146	31	M	Lung	adenocarcinoma, moderately differentiated, mucinous	T2N0M0	I B	6
147	62	M	Lung	adenocarcinoma, well differentiated, papillary	T2N0M0	B	6
148	63	М	Lung	adenocarcinoma, poorly differentiated	T2N0M0	I B	6
149	47	М	Lung	bronchioloalveolar carcinoma	T2N1M0	II B	6
150			Carbon				

TNM and Stage: AJCC Cancer Staging Manual (6th Edition)

	1	11	21	31	41	51	61	71	81	91	101	111	121	131	141
	2	20									10.00				
1	3	1													. 1
1	4	1		1				1					11		1
	5	8			45			5	8	95	83				145
staded area	6	12-1			3										
	7	10	1	1	-	H	-	-	1	-	-	-	1 1		- 22
	8	1	+		1	H		1		-					
	9	Se		· · · ·							20		3		- 22
	10	21						1		5.5	2		0-0		150