

■ 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

NBP2-30266 - Human Esophageal Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	Tissue type *
1	58	M	Esophagus	basaloid carcinoma	T3N0M0	II B	6
2	54	M	Esophagus	squamous cell carcinoma, moderately differentiated	T2N3M0	III C	6
3	57	M	Esophagus	squamous cell carcinoma, poorly differentiated	T3N2M0	III B	6
4	75	M	Esophagus	squamous cell carcinoma	T3N1M0	III A	6
5	46	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III A	6
6	58	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N2M0	III B	6
7	70	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N2M0	III B	6
8	66	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0	II B	6
9	53	M	Esophagus	small cell carcinoma	T2N2M0	III A	6
10	67	M	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	II A	6
11	60	F	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	I B	6
12	69	M	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	II A	6
13	62	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0	II B	6
14	57	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III A	6
15	62	M	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	II A	6
16	36	F	Esophagus	adenocarcinoma, well differentiated	T1bN0M0	I A	6
17	60	M	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	I B	6
18	50	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0	II A	6
19	60	F	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0	III A	6
20	57	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N2M0	III B	6
21	58	M	Esophagus	squamous cell carcinoma, well differentiated	T3N2M0	III B	6
22	74	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0	II B	6
23	56	F	Esophagus	squamous cell carcinoma, poorly differentiated	T3N2M0	III B	6
24	43	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III A	6
25	69	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N2M0	III B	6
26	58	M	Esophagus	squamous cell carcinoma, poorly differentiated	T3N1M0	III A	6
27	62	M	Esophagus	squamous cell carcinoma, well differentiated	T4aN0M0	III A	6
28	63	M	Esophagus	adenosquamous carcinoma	T3N1M0	III A	6
29	56	M	Esophagus	squamous cell carcinoma, well differentiated	T4aN2M0	III C	6
30	59	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0	III A	6
31	50	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III A	6
32	52	M	Esophagus	squamous cell carcinoma, well differentiated	T2N2M0	III A	6
33	58	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N3M0	III C	6
34	64	F	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0	III A	6
35	60	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0	III A	6
36	64	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M0	III A	6
37	52	M	Esophagus	squamous cell carcinoma, well differentiated	T3N2M0	III B	6
38	45	M	Esophagus	adenosquamous carcinoma	T3N0M0	I B	6
39	46	M	Esophagus	squamous cell carcinoma, well differentiated	T3N0M0	I B	6
40	66	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0	II B	6
41	69	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0	III A	6
42	60	M	Esophagus	squamous cell carcinoma, well differentiated	T3N2M0	III B	6
43	62	M	Esophagus	squamous cell carcinoma, well differentiated	T3N1M1	IV	6
44	59	M	Esophagus	squamous cell carcinoma, poorly differentiated	T3N1M0	III A	6
45	52	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0	II A	6
46	67	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0	III A	6
47	65	M	Esophagus	adenosquamous cell carcinoma	T4aN0M0	III A	6
48	51	M	Esophagus	adenosquamous cell carcinoma, moderately differentiated	T3N0M0	II B	6
49	47	M	Esophagus	squamous cell carcinoma, well differentiated	T3N2M0	III B	6
50	42	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N3M1	IV	6
51	66	M	Esophagus	squamous cell carcinoma, poorly differentiated	T3N1M0	III A	6
52	53	M	Esophagus	adenocarcinoma, moderately differentiated	T3N2M0	III B	6
53	59	M	Esophagus	squamous cell carcinoma, well differentiated	T3N2M0	III B	6
54	59	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0	III A	6
55	50	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N1M0	III A	6
56	42	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0	II A	6
57	69	M	Esophagus	adenocarcinoma, poorly differentiated	T3N3M0	III C	6
58	54	M	Esophagus	squamous cell carcinoma, moderately differentiated	T3N0M0	II A	6
59	70	M	Esophagus	squamous cell carcinoma, moderately differentiated	T2N1M0	II B	6
60	68	M	Esophagus	squamous cell carcinoma, well differentiated	T3N2M0	III B	6

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