

## ■ 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

## NBP2-30277 - Human Lung Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	LN*	pTNM	Stage	No. of NBP2-30277#	Tissue type**
1	56	M	Lung	squamous cell carcinoma, well differentiated	5/43	T2aN1M0	IIA	101	6
2	47	F	Lung	adenocarcinoma, moderately differentiated	0/30	T2aN0M0	IB	102	6
3	34	F	Lung	adenocarcinoma, well differentiated	18/30	T2bN2M0	IIIA	103	6
4	69	F	Lung	adenocarcinoma, well differentiated	0/15	T2aN0M0	IB	104	6
5	59	M	Lung	squamous cell carcinoma, moderately differentiated	1/15	T2bN1M0	IIB	105	6
6	62	M	Lung	squamous cell carcinoma, moderately differentiated	0/8	T2aN0M0	IB	106	6
7	42	M	Lung	bronchioloalveolar carcinoma, non-mucinous	7/17	T3N2M0	IIIA	107	6
8	64	M	Lung	adenocarcinoma, well differentiated	0/39	T2aN0M0	IB	108	6
9	64	M	Lung	squamous cell carcinoma, well differentiated	0/15	T2aN0M0	IB	109	6
10	68	M	Lung	squamous cell carcinoma, moderately differentiated	0/16	T3N0M0	IIB	110	6
11	54	M	Lung	carcinosarcoma	0/6	T3N0M0	IIB	111	6
12	73	M	Lung	large cell neuroendocrine carcinoma	0/7	T2aN0M0	IB	112	6
13	66	M	Lung	squamous cell carcinoma, well differentiated	2/20	T3N2M0	IIIA	113	6
14	59	F	Lung	large cell carcinoma	0/9	T2aN0M0	IB	114	6
15	59	M	Lung	squamous cell carcinoma, well differentiated	4/61	T2aN1M0	IIA	115	6
16	65	M	Lung	bronchioloalveolar carcinoma, non-mucinous	0/8	T2bN0M0	IIA	116	6
17	64	M	Lung	bronchioloalveolar carcinoma, mucinous	0/8	T2aN0M0	IB	117	6
18	61	M	Lung	squamous cell carcinoma, moderately differentiated	0/16	T2aN0M0	IB	118	6
19	69	F	Lung	squamous cell carcinoma, moderately differentiated	0/34	T2bN0M0	IIA	119	6
20	64	M	Lung	squamous cell carcinoma, well differentiated	0/17	T2aN0M0	IB	120	6
21	74	M	Lung	squamous cell carcinoma, well differentiated	1/46	T2bN1M0	IIB	121	6
22	53	F	Lung	squamous cell carcinoma, spindle cell	0/10	T1bN0M0	IA	122	6
23	61	M	Lung	squamous cell carcinoma, well differentiated	1/43	T2aN1M0	IIA	123	6
24	59	F	Lung	bronchioloalveolar carcinoma, mucinous	0/24	T2aN0M0	IB	124	6
25	58	M	Lung	large cell carcinoma	0/14	T1bN0M0	IA	125	6
26	77	M	Lung	squamous cell carcinoma, moderately differentiated	5/18	T2aN2M0	IIIA	126	6
27	53	F	Lung	bronchioloalveolar carcinoma, mucinous	0/9	T3N0M0	IIB	127	6
28	72	M	Lung	squamous cell carcinoma, poorly differentiated	0/18	T2bN0M0	IIA	128	6
29	63	M	Lung	squamous cell carcinoma, moderately differentiated	1/25	T2aN1M0	IIA	129	6
30	63	M	Lung	squamous cell carcinoma, moderately differentiated	0/22	T2bN0M0	IIA	130	6
31	59	M	Lung	squamous cell carcinoma, moderately differentiated	3/21	T2aN1M0	IIA	131	6
32	62	M	Lung	squamous cell carcinoma, poorly differentiated	0/50	T4N0M0	IIIA	132	6
33	66	M	Lung	large cell neuroendocrine carcinoma	0/38	T3N0M0	IIB	133	6
34	58	M	Lung	squamous cell carcinoma, moderately differentiated	2/18	T1bN1M0	IIA	134	6
35	55	F	Lung	squamous cell carcinoma, poorly differentiated	1/3	T2aN2M0	IIIA	135	6
36	63	M	Lung	squamous cell carcinoma, well differentiated	8/28	T2aN2M0	IIIA	136	6
37	81	M	Lung	squamous cell carcinoma, well differentiated	0/17	T2aN0M0	IB	137	6
38	54	M	Lung	squamous cell carcinoma, moderately differentiated	1/33	T2aN1M0	IIA	138	6
39	49	M	Lung	squamous cell carcinoma, well differentiated	0/19	T2aN0M0	IB	139	6
40	68	M	Lung	squamous cell carcinoma, poorly differentiated	0/22	T3N0M0	IIB	140	6
41	60	M	Lung	squamous cell carcinoma, poorly differentiated	0/22	T2aN0M0	IB	141	6
42	65	M	Lung	large cell carcinoma	0/33	T2aN0M0	IB	142	6
43	51	M	Lung	bronchioloalveolar carcinoma, mucinous	1/11	T1bN1M0	IIA	143	6
44	56	M	Lung	squamous cell carcinoma, well differentiated	4/32	T2aN2M0	IIIA	144	6
45	46	M	Lung	squamous cell carcinoma, moderately differentiated	1/25	T2aN1M0	IIA	145	6
46	71	M	Lung	squamous cell carcinoma, moderately differentiated	0/21	T2aN0M0	IB	146	6
47	69	M	Lung	squamous cell carcinoma, poorly differentiated	0/11	T2aN0M0	IB	147	6
48	58	M	Lung	squamous cell carcinoma, well differentiated	0/24	T2bN0M0	IIA	148	6
49	62	M	Lung	bronchioloalveolar carcinoma, non-mucinous	2/24	T1aN1M0	IIA	149	6
50	71	M	Lung	squamous cell carcinoma, well differentiated	5/40	T2bN1M0	IIB	150	6
51	65	M	Lung	squamous cell carcinoma, moderately differentiated	0/28	T3N0M0	IIB	151	6
52	67	M	Lung	squamous cell carcinoma, poorly differentiated	0/12	T2aN0M0	IB	152	6
53	64	M	Lung	squamous cell carcinoma, poorly differentiated	0/14	T3N0M0	IIB	153	6
54	63	M	Lung	carcinosarcoma	0/14	T2aN0M0	IB	154	6
55	69	M	Lung	squamous cell carcinoma, moderately differentiated	0/20	T2aN0M0	IB	155	6
56	33	F	Lung	adenocarcinoma, moderately differentiated	0/6	T2aN0M0	IB	156	6
57	60	F	Lung	adenocarcinoma, well differentiated	22/22	T3N2M0	IIIA	157	6
58	68	M	Lung	adenocarcinoma, moderately differentiated	0/12	T2aN0M0	IB	158	6
59	41	F	Lung	adenocarcinoma, moderately differentiated	2/15	T3N2M0	IIIA	159	6
60	.	.	Carbon	.	.	.	.	.	.

\*positive lymph nodes / examined lymph nodes

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

# : The normal tissue in NBP2-30223 of corresponding number is from the identical patient.