

## ■ 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

## NBP2-30231 - Human Uterine Cervix Tissue MicroArray (Cancer metastasis)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	Tissue type*
1	70	F	Cervix	squamous cell carcinoma, microinvasive	T1a1NOM0	IA1	6
2	46	F	Cervix	squamous cell carcinoma, microinvasive	T1a1NOM0	IA1	6
3	42	F	Cervix	squamous cell carcinoma, microinvasive	T1a1NOM0	IA1	6
4	57	F	Cervix	squamous cell carcinoma in situ	TisNOM0	0	6
5	33	F	Cervix	squamous cell carcinoma, microinvasive	T1a1NOM0	IA1	6
6	35	F	Cervix	squamous cell carcinoma in situ	TisNOM0	0	6
7	48	F	Cervix	squamous cell carcinoma, microinvasive	T1a1NOM0	IA1	6
8	39	F	Cervix	squamous cell carcinoma, microinvasive	T1a1NOM0	IA1	6
9	58	F	Cervix	squamous cell carcinoma, microinvasive	T1a1NOM0	IA1	6
10	53	F	Cervix	squamous cell carcinoma in situ	TisNOM0	0	6
11	53	F	Cervix	squamous cell carcinoma	T1b2NOM0	IB2	6
12	53	F	Cervix	squamous cell carcinoma	T2bN1M0	III B	6
13	50	F	Cervix	adenocarcinoma	T1b2N1M0	III B	6
14	45	F	Cervix	squamous cell carcinoma	T1b2N1M0	III B	6
15	49	F	Cervix	squamous cell carcinoma	T1b1N1M0	III B	6
16	35	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
17	57	F	Cervix	squamous cell carcinoma	T2a1NOM0	IIA1	6
18	39	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
19	66	F	Cervix	squamous cell carcinoma	T2a1N1M0	III B	6
20	55	F	Cervix	squamous cell carcinoma	T1b2NOM0	IB2	6
21	64	F	Cervix	squamous cell carcinoma	T1b2N1M0	III B	6
22	50	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
23	48	F	Cervix	squamous cell carcinoma	T1b1N1M0	III B	6
24	51	F	Cervix	squamous cell carcinoma	T1b2N1M0	III B	6
25	54	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
26	39	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
27	58	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
28	57	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
29	55	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
30	35	F	Cervix	squamous cell carcinoma	T1b1N1M0	III B	6
31	62	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
32	64	F	Cervix	squamous cell carcinoma	T1b2N1M0	III B	6
33	60	F	Cervix	adenosquamous carcinoma	T1b1NOM0	IB1	6
34	43	F	Cervix	squamous cell carcinoma	T1b1N1M0	III B	6
35	68	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
36	45	F	Cervix	squamous cell carcinoma	T1b2NOM0	IB2	6
37	40	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
38	58	F	Cervix	adenosquamous carcinoma	T1b1NOM0	IB1	6
39	67	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
40	48	F	Cervix	squamous cell carcinoma	T1b1N1M0	III B	6
41	62	F	Cervix	squamous cell carcinoma	T1b2N1M0	III B	6
42	32	F	Cervix	squamous cell carcinoma	T1b1N1M0	III B	6
43	55	F	Cervix	squamous cell carcinoma	T2a1N1M0	III B	6
44	61	F	Cervix	squamous cell carcinoma	T2bN1M0	III B	6
45	34	F	Cervix	adenocarcinoma	T1b2N1M0	III B	6
46	53	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
47	45	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
48	65	F	Cervix	squamous cell carcinoma	T1b1NOM0	IB1	6
49	47	F	Cervix	squamous cell carcinoma	T1b1N1M0	III B	6
50	59	F	Cervix	adenosquamous carcinoma	T1b2N1M0	III B	6
51	62	F	Lymph node	metastatic carcinoma of No.41	.	.	6
52	32	F	Lymph node	metastatic carcinoma of No.42	.	.	6
53	55	F	Lymph node	metastatic carcinoma of No.43	.	.	6
54	61	F	Lymph node	metastatic carcinoma of No.44	.	.	6
55	34	F	Lymph node	metastatic carcinoma of No.45	.	.	6
56	53	F	Cervix	normal of No.46	.	.	3
57	45	F	Cervix	normal of No.47	.	.	3
58	65	F	Cervix	normal of No.48	.	.	3
59	47	F	Cervix	normal of No.49	.	.	3
60	.	.	Carbon	.	.	.	.

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