

■ 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

NBP2-30218 - Human Endometrium Tissue MicroArray (Normal)

No.	Age	Sex	Organ	Diagnosis	Associated lesion	Tissue type*
1	35	F	Endometrium	proliferative phase	squamous cell carcinoma of cervix	3
2	39	F	Endometrium	proliferative phase	leiomyoma	1
3	40	F	Endometrium	proliferative phase	squamous cell carcinoma in situ of cervix	3
4	40	F	Endometrium	secretory phase	leiomyoma	1
5	33	F	Endometrium	secretory phase	leiomyoma	1
6	45	F	Endometrium	secretory phase	leiomyoma	1
7	40	F	Endometrium	secretory phase	leiomyoma	1
8	39	F	Endometrium	secretory phase	leiomyoma	1
9	38	F	Endometrium	proliferative phase	endometriosis, ovary	1
10	42	F	Endometrium	secretory phase	leiomyoma	1
11	44	F	Endometrium	secretory phase	leiomyoma	1
12	45	F	Endometrium	secretory phase	leiomyoma	1
13	37	F	Endometrium	secretory phase	leiomyoma	1
14	38	F	Endometrium	proliferative phase	leiomyoma	1
15	42	F	Endometrium	secretory phase	squamous cell carcinoma of cervix	3
16	38	F	Endometrium	proliferative phase	leiomyoma	1
17	37	F	Endometrium	proliferative phase	leiomyoma	1
18	41	F	Endometrium	proliferative phase	leiomyoma	1
19	34	F	Endometrium	secretory phase	squamous cell carcinoma in situ of cervix	3
20	43	F	Endometrium	secretory phase	leiomyoma	1
21	40	F	Endometrium	proliferative phase	leiomyoma	1
22	42	F	Endometrium	secretory phase	leiomyoma	1
23	41	F	Endometrium	proliferative phase	moderate dysplasia of cervix	1
24	45	F	Endometrium	proliferative phase	leiomyoma	1
25	44	F	Endometrium	proliferative phase	leiomyoma	1
26	30	F	Endometrium	secretory phase	squamous cell carcinoma of cervix	3
27	39	F	Endometrium	proliferative phase	leiomyoma	1
28	43	F	Endometrium	secretory phase	leiomyoma	1
29	36	F	Endometrium	proliferative phase	leiomyoma	1
30	44	F	Endometrium	secretory phase	leiomyoma	1
31	42	F	Endometrium	proliferative phase	leiomyoma	1
32	30	F	Endometrium	secretory phase	leiomyoma	1
33	41	F	Endometrium	secretory phase	leiomyoma	1
34	33	F	Endometrium	secretory phase	leiomyoma	1
35	37	F	Endometrium	secretory phase	normal	1
36	29	F	Endometrium	secretory phase	squamous cell carcinoma in situ of cervix	3
37	42	F	Endometrium	secretory phase	leiomyoma	1
38	37	F	Endometrium	proliferative phase	squamous cell carcinoma in situ of cervix	3
39	44	F	Endometrium	proliferative phase	leiomyoma	1
40	35	F	Endometrium	proliferative phase	adenomyoma	1
41	32	F	Endometrium	secretory phase	leiomyoma	1
42	36	F	Endometrium	secretory phase	cellular leiomyoma	1
43	37	F	Endometrium	secretory phase	leiomyoma	1
44	35	F	Endometrium	secretory phase	leiomyoma	1
45	37	F	Endometrium	secretory phase	squamous cell carcinoma in situ of cervix	3
46	42	F	Endometrium	secretory phase	leiomyoma	1
47	44	F	Endometrium	proliferative phase	leiomyoma	1
48	40	F	Endometrium	proliferative phase	cellular leiomyoma	1
49	35	F	Endometrium	proliferative phase	squamous cell carcinoma of cervix	3
50	36	F	Endometrium	proliferative phase	squamous cell carcinoma of cervix	3
51	35	F	Endometrium	secretory phase	squamous cell carcinoma of cervix	3
52	43	F	Endometrium	secretory phase	squamous cell carcinoma of cervix	3
53	42	F	Endometrium	proliferative phase	leiomyoma	1
54	35	F	Endometrium	proliferative phase	squamous cell carcinoma of cervix	3
55	43	F	Endometrium	secretory phase	leiomyoma	1
56	42	F	Endometrium	secretory phase	squamous cell carcinoma of cervix	3
57	33	F	Endometrium	proliferative phase	leiomyoma	1
58	35	F	Endometrium	proliferative phase	adenomyosis	1
59	45	F	Endometrium	secretory phase	leiomyoma	1
60	.	.	Carbon	.	.	.