

■ 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

NBP2-30290 - Human Ovary Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	Tissue type*
1	65	F	Ovary	papillary serous tumor of borderline malignancy	T1aNOM0	IA	5
2	26	F	Ovary	mucinous cystadenoma of borderline malignancy	T1aNOM0	IA	5
3	35	F	Ovary	papillary serous tumor of borderline malignancy	T1aNOM0	IA	5
4	65	F	Ovary	serous cystadenocarcinoma	T2bNOM0	IIB	6
5	48	F	Ovary	endometrioid carcinoma	T2bNOM0	IIB	6
6	44	F	Ovary	granulosa cell tumor	T1aNOM0	IA	6
7	49	F	Ovary	endometrioid carcinoma	T2bNOM0	IIB	6
8	67	F	Ovary	mucinous cystadenoma of borderline malignancy	T2bNOM0	IIB	5
9	75	F	Ovary	serous cystadenofibroma	T1aNOM0	IA	4
10	50	F	Ovary	papillary serous cystadenocarcinoma , poorly differentiated	T2cNOM0	IIC	6
11	44	F	Ovary	papillary serous tumor of borderline malignancy	T1aNOM0	IA	5
12	50	F	Ovary	papillary serous cystadenocarcinoma	T3cNOM1	IV	6
13	47	F	Ovary	endometrioid carcinoma	T3cNOM1	IV	6
14	38	F	Ovary	mixed Brenner tumor and mucinous cystadenoma	T1aNOM0	IA	4
15	58	F	Ovary	papillary serous cystadenocarcinoma	T2bNOM0	IIB	6
16	70	F	Ovary	papillary serous cystadenocarcinoma	T3bNOM0	IIIB	6
17	65	F	Ovary	struma ovarii	T1aNOM0	IA	4
18	79	F	Ovary	endometrioid adenoacanthofibroma	T1aNOM0	IA	4
19	57	F	Ovary	serous surface papillary carcinoma	T2bNOM0	IIB	6
20	43	F	Ovary	papillary serous cystadenocarcinoma	T3bNOM0	IIIB	6
21	68	F	Ovary	undifferentiated carcinoma	T1bNOM0	IB	6
22	50	F	Ovary	clear cell carcinoma	T1aNOM0	IA	6
23	23	F	Ovary	dysgerminoma	T2bNOM0	IIB	6
24	51	F	Ovary	common epithelial carcinoma , poorly differentiated	T2bNOM0	IIB	6
25	70	F	Ovary	malignant muellerian mixed tumor	T1aNOM0	IA	6
26	21	F	Ovary	dysgerminoma	T1aNOM0	IA	6
27	43	F	Ovary	papillary serous cystadenocarcinoma	T2bN1bM0	IIIA1	6
28	61	F	Ovary	papillary serous cystadenocarcinoma	T3bNOM0	IIIB	6
29	54	F	Ovary	sertoli-leydig cell tumor	T1bNOM0	IB	6
30	69	F	Ovary	granulosa cell tumor	T1aNOM0	IA	6
31	38	F	Ovary	clear cell carcinoma	T1aNOM1b	IVB	6
32	18	F	Ovary	yolk sac tumor	T1aNOM0	IA	6
33	31	F	Ovary	dysgerminoma	T2cNOM0	IIC	6
34	51	F	Ovary	papillary serous cystadenocarcinoma	T2bN1M0	IIIC	6
35	61	F	Ovary	granulosa cell tumor	T1aNOM0	IA	6
36	24	F	Ovary	sertoli-leydig cell tumor	T1aNOM0	IA	6
37	15	F	Ovary	yolk sac tumor	T1aNOM0	IA	6
38	41	F	Ovary	clear cell carcinoma	T2bNOM0	IIB	6
39	56	F	Ovary	malignant muellerian mixed tumor	T1aNOM0	IA	6
40	41	F	Ovary	clear cell carcinoma	T1aNOM0	IA	6
41	27	F	Ovary	embryonal carcinoma	T2aNOM0	IIA	6
42	55	F	Ovary	granulosa cell tumor	T1aNOM0	IA	5
43	23	F	Ovary	dysgerminoma	T1aNOM0	IA	6
44	64	F	Ovary	papillary serous cystadenocarcinoma	T2cN1M0	IIIC	6
45	43	F	Ovary	metastatic adenocarcinoma (most likely from breast)	M1	IV	6
46	48	F	Ovary	malignant lymphoma	.	II	6
47	50	F	Ovary	adenocarcinoma from salpinx	T2bNOM0	IIB	6
48	16	F	Ovary	mixed germ cell tumor	T1bN1aM0	IIIA1	6
49	46	F	Ovary	metastatic undifferentiated carcinoma	M1	IV	6
50	37	F	Ovary	metastatic signet ring carcinoma from stomach	M1	IV	6
51	55	F	Ovary	serous cystadenocarcinoma, moderately differentiated	T3bNOM0	IIIB	6
52	18	F	Ovary	dysgermonima	T1aNOM0	IA	6
53	45	F	Ovary	serous cystadenocarcinoma, moderately differentiated	T2cNOM0	IIC	6
54	37	F	Ovary	endometrioid carcinoma	T2bNOM0	IIB	6
55	14	F	Ovary	yolk sac tumor	T1cNOM0	IC	6
56	65	F	Ovary	adenocarcinoma, poorly differentiated	T3cNOM0	IIIC	6
57	39	F	Ovary	clear cell carcinoma	T2cN1M0	IIIC	6
58	40	F	Ovary	papillary serous cystadenocarcinoma	T1aNOM0	IA	6
59	58	F	Ovary	clear cell carcinoma	T1aNOM0	IA	6
60	.	.	Carbon

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