

■ 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

NBP2-30317 - Human Thyroid Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	Tissue type*
1	38	F	Thyroid	papillary carcinoma	T3N1aM0	I	6
2	43	F	Thyroid	papillary carcinoma	T3N0M0	I	6
3	31	F	Thyroid	papillary carcinoma	T3N1aM0	I	6
4	35	F	Thyroid	papillary carcinoma	T3N0M0	I	6
5	41	F	Thyroid	papillary carcinoma	T3N0M0	I	6
6	57	F	Thyroid	papillary carcinoma	T3N0M0	III	6
7	52	F	Thyroid	papillary carcinoma	T2N0M0	II	6
8	45	M	Thyroid	papillary carcinoma	T3N0M0	III	6
9	40	F	Thyroid	papillary carcinoma	T3N0M0	I	6
10	60	F	Thyroid	papillary carcinoma	T3N1aM0	III	6
11	48	F	Thyroid	papillary carcinoma	T3N0M0	III	6
12	29	F	Thyroid	papillary carcinoma	T2N1aM0	I	6
13	72	M	Thyroid	papillary carcinoma	T3N1bM0	IVA	6
14	24	M	Thyroid	papillary carcinoma	T3N1aM0	I	6
15	37	F	Thyroid	follicular carcinoma	T2N0M0	I	6
16	56	F	Thyroid	papillary carcinoma	T3N1bM0	IVA	6
17	35	M	Thyroid	follicular carcinoma	T3N0M0	I	6
18	55	F	Thyroid	papillary carcinoma	T3N0M0	III	6
19	35	F	Thyroid	papillary carcinoma	T3N0M0	I	6
20	54	M	Thyroid	papillary carcinoma	T3N1aM0	III	6
21	28	F	Thyroid	papillary carcinoma	T3N1bM0	I	6
22	45	F	Thyroid	papillary carcinoma	T3N1aM0	III	6
23	46	M	Thyroid	papillary carcinoma	T3N0M0	III	6
24	36	F	Thyroid	papillary carcinoma	T3N1aM0	I	6
25	67	F	Thyroid	papillary carcinoma	T3N1aM0	III	6
26	22	F	Thyroid	follicular carcinoma	T2N0M0	I	6
27	45	M	Thyroid	papillary carcinoma	T3N1aM0	III	6
28	33	F	Thyroid	follicular carcinoma	T2N0M0	I	6
29	36	F	Thyroid	papillary carcinoma	T2N0M0	I	6
30	21	F	Thyroid	papillary carcinoma	T2N0M0	I	6
31	22	F	Thyroid	papillary carcinoma	T3N1aM0	I	6
32	18	F	Thyroid	papillary carcinoma	T3N1aM0	I	6
33	17	F	Thyroid	papillary carcinoma	T2N0M0	I	6
34	51	F	Thyroid	papillary carcinoma	T3N1aM0	III	6
35	50	F	Lymph node	papillary carcinoma, recurrent	rTON1bM0	IVA	6
36	57	F	Thyroid	follicular carcinoma	T2N0M0	II	6
37	38	F	Thyroid	papillary carcinoma	T3N1bM0	I	6
38	20	F	Thyroid	papillary carcinoma	T3N1aM0	I	6
39	27	F	Thyroid	follicular carcinoma	T2N0M0	I	6
40	65	M	Thyroid	follicular carcinoma	T2N0M0	II	6
41	40	F	Thyroid	papillary carcinoma	T3N0M0	I	6
42	69	F	Thyroid	papillary carcinoma	T3N1aM0	III	6
43	50	F	Thyroid	papillary carcinoma	T3N0M0	III	6
44	51	M	Thyroid	follicular carcinoma	T2N0M0	II	6
45	53	M	Thyroid	anaplastic carcinoma	T4aN0M0	IVA	6
46	47	M	Thyroid	papillary carcinoma	T3N1bM0	IVA	6
47	63	F	Thyroid	papillary carcinoma	T2N0M0	II	6
48	58	M	Thyroid	papillary carcinoma	T3N0M0	III	6
49	75	F	Thyroid	papillary carcinoma	T3N0M0	III	6
50	15	F	Thyroid	papillary carcinoma	T3N0M0	I	6
51	45	M	Thyroid	benign thyroid tissue (same patient of #8)	.	.	3
52	31	F	Thyroid	benign thyroid tissue	.	.	3
53	38	F	Thyroid	benign thyroid tissue	.	.	3
54	50	F	Thyroid	benign thyroid tissue (same patient of #43)	.	.	3
55	52	M	Thyroid	benign thyroid tissue	.	.	3
56	78	F	Thyroid	benign thyroid tissue	.	.	3
57	32	F	Thyroid	benign thyroid tissue	.	.	3
58	15	F	Thyroid	benign thyroid tissue (same patient of #50)	.	.	3
59	37	F	Thyroid	benign thyroid tissue	.	.	3
60	.	.	Carbon

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