

■ 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

NBP2-30229 - Human Skin Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Tissue type*
1	80	M	Skin, leg	invasive squamous cell carcinoma, well differentiated	T4N0M0	6
2	64	M	Skin, lower leg	squamous cell carcinoma	T4N0M0	6
3	62	M	Skin, neck	squamous cell carcinoma, well differentiated	T2N0M0	6
4	58	F	Skin, foot	squamous cell carcinoma, well differentiated	T2N0M0	6
5	55	M	Skin, neck	squamous cell carcinoma, moderately differentiated	T4N0M0	6
6	60	F	Skin, face	squamous cell carcinoma, moderately differentiated	T2N0M0	6
7	58	M	Skin, shin	squamous cell carcinoma, well differentiated	T4N0M0	6
8	54	M	Skin, lower leg	squamous cell carcinoma, well differentiated	T4N0M0	6
9	59	M	Skin, sacral area	squamous cell carcinoma, well differentiated	T1N0M0	6
10	67	M	Skin, neck	squamous cell carcinoma, well differentiated, recurrent	rT2N0M0	6
11	64	F	Skin, ear	basal cell carcinoma	T2N0M0	6
12	46	M	Skin, inguinal area	squamous cell carcinoma	T2N1M0	6
13	54	F	Skin, buttock	squamous cell carcinoma, moderately differentiated	T2N0M0	6
14	39	F	Skin, forearm	squamous cell carcinoma, well differentiated	T2N0M0	6
15	65	M	Skin, cheek	squamous cell carcinoma, well differentiated	T2N0M0	6
16	65	M	Skin, perianal	squamous cell carcinoma, well differentiated	T2N0M0	6
17	46	M	Skin, lower leg	invasive squamous cell carcinoma, well differentiated	T4N0M0	6
18	68	F	Skin, thumb	squamous cell carcinoma	T1N0M0	6
19	58	M	Skin, lip	squamous cell carcinoma, well differentiated	T3N1M0	6
20	71	M	Skin, lower leg	squamous cell carcinoma, well differentiated	T4N0M0	6
21	40	M	Skin, lower leg	squamous cell carcinoma, well differentiated	T2N0M0	6
22	65	M	Skin, left malleolus	squamous cell carcinoma	T2N0M0	6
23	76	F	Skin, scalp	squamous cell carcinoma, well differentiated	T2N0M0	6
24	75	F	Skin, scalp	squamous cell carcinoma, well differentiated	T4N0M0	6
25	66	M	Skin, lower lip	squamous cell carcinoma, moderately differentiated	T1N2bM0	6
26	95	F	Skin, thigh	basal cell carcinoma, mixed type	T2N0M0	6
27	55	M	Skin, foot	verrucous carcinoma	T2N0M0	6
28	72	M	Skin, neck	squamous cell carcinoma, well differentiated	T2N0M0	6
29	36	F	Vulva	microinvasive squamous cell carcinoma, well differentiated	T1bN0M0	6
30	45	M	Skin, planta	squamous cell carcinoma, well differentiated	T2N0M0	6
31	71	F	Skin, scalp	basal cell carcinoma with adnexal differentiation	T2N0M0	6
32	85	F	Skin, cheek	squamous cell carcinoma, well differentiated	T1N0M0	6
33	67	M	Skin, neck	squamous cell carcinoma, well differentiated	T1N1M0	6
34	42	M	Skin, thigh	squamous cell carcinoma, recurrent	rT2N1M1	6
35	63	M	Skin, chest wall	squamous cell carcinoma, moderately differentiated	T2N1M1	6
36	56	M	Skin, ear	basal cell carcinoma	T1N0M0	6
37	64	M	Skin, arm	squamous cell carcinoma, moderately differentiated	T2N0M0	6
38	74	F	Skin, face	basal cell carcinoma	T1N0M0	6
39	42	M	Tongue	squamous cell carcinoma, well differentiated	T3N1M0	6
40	71	M	Skin, flank	Bowen's disease	Tis	6
41	65	M	Skin	normal	.	3
42	55	M	Skin	normal	.	3
43	64	M	Skin	normal of No. 37	.	3
44	76	F	Skin	normal of No. 23	.	3
45	65	M	Skin	normal	.	3
46	55	M	Skin	normal	.	3
47	80	F	Skin	normal	.	3
48	58	F	Skin	normal	.	3
49	65	F	Skin	normal	.	3
50	.	.	Carbon	.	.	.

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