

■ 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

NBP2-30215 - Human Multi Tissue MicroArray (Normal Adjacent)

No.	Age	Sex	Organ	Diagnosis	No. of NBP2-30262*	Tissue type**
101	56	M	Stomach	.	1	3
102	52	M	Stomach	.	2	3
103	39	M	Stomach	.	3	3
104	44	M	Stomach	.	4	3
105	65	F	Stomach	.	5	3
106	52	M	Stomach	.	6	3
107	66	M	Stomach	.	7	3
108	61	M	Stomach	.	8	3
109	39	F	Stomach	.	9	3
110	60	F	Stomach	.	10	3
111	64	M	Esophagus	.	11	3
112	60	M	Esophagus	.	12	3
113	73	M	Esophagus	.	13	3
114	52	F	Esophagus	.	14	3
115	77	M	Esophagus	.	15	3
116	63	M	Esophagus	.	16	3
117	47	M	Esophagus	.	17	3
118	58	M	Esophagus	.	18	3
119	60	F	Esophagus	.	19	3
120	69	M	Esophagus	.	20	3
121	69	M	Lung	.	21	3
122	71	M	Lung	.	22	3
123	65	M	Lung	.	23	3
124	62	M	Lung	.	24	3
125	69	M	Lung	.	25	3
126	58	M	Lung	.	26	3
127	58	M	Lung	.	27	3
128	61	M	Lung	.	28	3
129	52	M	Lung	.	29	3
130	56	M	Lung	.	30	3
131	66	M	Sigmoid colon	.	31	3
132	50	M	Rectum	.	32	3
133	40	M	Cecum	.	33	3
134	45	M	Ascending colon	.	34	3
135	43	F	Rectum	.	35	3
136	44	F	Ascending colon	.	36	3
137	52	F	Rectum	.	37	3
138	49	M	Ascending colon	.	38	3
139	61	F	Rectum	.	39	3
140	52	F	Rectum	.	40	3
141	54	F	Thyroid	.	.	3
142	20	F	Thyroid	.	42	3
143	15	F	Thyroid	.	43	3
144	60	F	Thyroid	.	44	3
145	47	F	Thyroid	.	45	3
146	48	F	Thyroid	.	46	3
147	43	F	Thyroid	.	47	3
148	56	F	Thyroid	.	48	3
149	55	F	Thyroid	.	49	3
150	34	F	Thyroid	.	50	3
151	57	F	Kidney	.	51	3
152	50	F	Kidney	.	52	3
153	27	M	Kidney	.	53	3
154	56	M	Kidney	.	54	3
155	47	M	Kidney	.	55	3
156	63	M	Kidney	.	56	3
157	53	M	Kidney	.	57	3
158	62	M	Kidney	.	58	3
159	65	F	Kidney	.	.	3
160	.	.	Carbon	.	.	.

*: The cancer tissue in NBP2-30262 of corresponding number is from the identical patient.