

■ 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
	31	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

NBP2-30216 - Human Multi Tissue MicroArray (Normal Adjacent)

No.	Age	Sex	Organ	Diagnosis	No. of NBP2-30263*	Tissue type**
101	30	F	Breast	.	.	3
102	43	F	Breast	.	2	3
103	58	F	Breast	.	3	3
104	46	F	Breast	.	4	3
105	71	F	Breast	.	5	3
106	37	F	Breast	.	6	3
107	31	F	Breast	.	.	3
108	30	F	Breast	.	8	3
109	47	F	Breast	.	9	3
110	49	F	Breast	.	.	3
111	32	M	Liver	chronic active hepatitis	11	3
112	57	M	Liver	cirrhosis	12	3
113	59	F	Liver	chronic active hepatitis	.	3
114	26	M	Liver	biliary cirrhosis	.	3
115	62	M	Liver	non-specific reactive change	.	3
116	64	M	Liver	chronic active hepatitis	.	3
117	61	M	Liver	non-specific reactive change	17	3
118	60	M	Liver	cirrhosis	.	3
119	50	M	Liver	chronic active hepatitis	19	3
120	62	M	Liver	cirrhosis	.	3
121	53	M	Skeletal muscle	.	21	2
122	52	F	Urinary bladder, mucosa	.	.	3
123	60	M	Lymph node	.	23	2
124	65	M	Urinary bladder, mucosa	.	24	3
125	74	M	Prostate	.	25	2
126	63	F	Urinary bladder, mucosa	.	26	3
127	55	M	Urinary bladder, mucosa	.	.	3
128	46	M	Urinary bladder, mucosa	.	28	3
129	52	M	Urinary bladder, mucosa	.	.	3
130	63	M	Prostate	.	30	2
131	46	F	Ovary	.	.	3
132	50	F	Ovary	.	.	3
133	60	F	Endometrium	.	33	2
134	41	F	Ovary	.	.	3
135	64	F	Myometrium	.	35	2
136	46	F	Ovary	.	.	3
137	35	F	Ovary	.	.	2
138	34	F	Ovary	.	.	3
139	37	F	Ovary	.	.	3
140	37	F	Ovary	.	.	3
141	66	M	Duodenum, mucosa	.	41	2
142	69	M	Duodenum, mucosa	.	.	2
143	68	M	Duodenum, mucosa	.	43	2
144	45	M	Duodenum, mucosa	.	44	2
145	66	M	Pancreas	.	.	3
146	29	F	Pancreas	.	.	3
147	69	M	Spleen	.	47	2
148	27	F	Pancreas	.	.	3
149	63	M	Pancreas	.	.	3
150	40	M	Pancreas	chronic pancreatitis	50	3
151	72	M	Prostate	.	.	3
152	76	M	Prostate	.	.	3
153	74	M	Prostate	.	.	3
154	74	M	Prostate	.	.	3
155	80	M	Prostate	.	.	3
156	68	M	Prostate	.	.	3
157	65	M	Prostate	.	.	3
158	74	M	Prostate	.	.	3
159	54	M	Prostate	.	59	3
160	.	.	Carbon	.	.	.

* : The cancer tissue in NBP2-30263 array of corresponding number is from the identical patient.