

■ 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

NBP2-30230 - Human Stomach Tissue MicroArray (Normal)

No.	Age	Sex	Organ	Diagnosis	No. of NBP2-30308#	Tissue type*
101	40	M	Stomach	.	1	3
102	43	F	Stomach	.	2	3
103	61	F	Stomach	.	3	3
104	45	F	Stomach	.	4	3
105	67	M	Stomach	.	5	3
106	53	F	Stomach	.	6	3
107	67	M	Stomach	.	7	3
108	72	M	Stomach	.	8	3
109	61	M	Stomach	.	9	3
110	65	M	Stomach	.	10	3
111	70	M	Stomach	.	11	3
112	62	M	Stomach	.	12	3
113	62	F	Stomach	.	11	3
114	52	M	Stomach	.	14	3
115	60	M	Stomach	.	15	3
116	66	M	Stomach	.	16	3
117	41	M	Stomach	.	17	3
118	60	F	Stomach	.	18	3
119	59	M	Stomach	.	19	3
120	62	M	Stomach	.	20	3
121	49	M	Stomach	.	21	3
122	64	M	Stomach	.	22	3
123	55	M	Stomach	.	23	3
124	56	F	Stomach	.	24	3
125	48	M	Stomach	.	25	3
126	65	F	Stomach	.	26	3
127	63	M	Stomach	.	27	3
128	60	M	Stomach	.	28	3
129	42	M	Stomach	.	29	3
130	55	M	Stomach	.	30	3
131	58	M	Stomach	.	31	3
132	72	F	Stomach	.	.	3
133	53	M	Stomach	.	33	3
134	46	M	Stomach	.	34	3
135	64	M	Stomach	.	35	3
136	65	M	Stomach	.	36	3
137	45	F	Stomach	.	37	3
138	58	F	Stomach	.	38	3
139	68	F	Stomach	.	39	3
140	66	M	Stomach	.	40	3
141	63	M	Stomach	.	41	3
142	52	F	Stomach	.	42	3
143	35	F	Stomach	.	43	3
144	44	F	Stomach	.	44	3
145	55	M	Stomach	.	45	3
146	40	M	Stomach	.	46	3
147	66	M	Stomach	.	47	3
148	65	F	Stomach	.	48	3
149	62	F	Stomach	.	49	3
150	60	F	Stomach	.	.	3
151	70	F	Stomach	.	51	3
152	52	M	Stomach	.	52	3
153	40	M	Stomach	.	53	3
154	34	F	Stomach	.	54	3
155	60	M	Stomach	.	55	3
156	68	M	Stomach	.	56	3
157	52	M	Stomach	.	57	3
158	62	M	Stomach	.	58	3
159	60	M	Stomach	.	59	3
160	.	.	Carbon	.	.	.

Tissue type*: see General information sheet

#: The cancer tissue in NBP2-30308 array of this column is from the identical patient.