

Data Sheet

Product Name: Frozen Tissue Panel

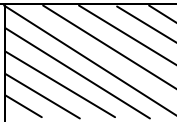
Catalog No.: **NBP2-78104**

Lot No.: **B806040**

Species: Human Mouse Rat ☒ Monkey (Rh) Guinea Pig Porcine
Bovine Hamster Dog Monkey (Cy) Rabbit Plant

Tissue Type: ☒ Normal Adult Fetal Tumor Disease Cell line

Product Image:

1. ☐	2. ☐	3. ☐	4. ☐	
5. ☐	6. ☐	7. ☐	8. ☐	

Donor Information

Item	Tissue Name	Sex	Age (Years)	Clinical Diagnosis
1	Heart	M	4	Normal
2	Brain	M	4.5	Normal
3	Kidney	M	2.5	Normal
4	Liver	M	2.5	Normal
5	Lung	M	2.5	Normal
6	Spleen	F	4.5	Normal
7	Small Intestine	F	4.5	Normal
8	Skeletal Muscle	M	2.5	Normal

Components

- 5 slides per package
- Certificate of Analysis

FOR IN VITRO RESEARCH USE ONLY

APPROVED BY: 