



DISCOVERY

L I F E S C I E N C E S

Science at your Service.TM

Recommended Handling and Storage for Biological Specimens

All Biological samples should be handling following Universal Precautions set forth in 29 CFR 1910.1030, Occupational Safety and Health Standards.

Ensure Proper Personal Protective Equipment (PPE) is worn at all times by all persons who come in contact with biological samples.

Cell Products (PBMCs, BMMCs, DTCs, Flash Frozen Tissue):

- For best results, store cryopreserved cell products in a liquid nitrogen freezer ($\leq -135^{\circ}\text{C}$) until ready to use. Follow Discovery's Thawing Protocol for best cell recovery results.

Non-cell Biologic Products (Plasma, Serum, Buffy Coat):

- For best results, store non-cell biological specimens in an Ultra-Low freezer ($\leq -80^{\circ}\text{C}$) until ready to use.

Formalin-Fixed Paraffin Embedded (FFPE) Tissue, Slides and Curls:

- Paraffin blocks can be stored at temperatures close to room temperature (15 to 25°C). Blocks should be stored out of direct sunlight and not in close proximity to an open flame or direct heat source.
- Slides cut from paraffin blocks can lose antigenicity when stored over long periods of time. This process can be slowed down by storing unused slides at 4°C . Loss of antigenicity can lead to suboptimal results, particularly when using these slides for Immunohistochemistry (IHC).
- Discovery Life Sciences recommends using slides within one calendar month from the date of sectioning.