

Cryo-EM Grid Storage and Shipping Puck Systems



Get informed about
[Pucks & Accessories ↗](#)



Get informed about
[Storage & Shipping ↗](#)



MiTeGen Cryo-EM Grid Storage and Shipping Puck Systems

allow you organize, store, track, and ship your samples on grids for single particle analysis and other Cryogenic Electron Microscopy (Cryo-EM) applications with ease and efficiency.

How it works - Your samples on grids are placed in grid boxes. The grid boxes are placed in a storage puck. The storage puck is placed into either a storage or shipping cane (rack). Then the cane is placed in either a storage dewar or dryshipper. You are now ready to store or ship your samples.

Tracking and organization - There are several options for tracking and keeping your samples organized. These include barcoding and/or serialization of the pucks, puck color selection, barcode scanning as well as log sheets.

Several system sizes - We offer system ranging from the basic system for those just getting started in cryo-EM to systems for facilities processing large number of samples.

We offer several cryo-EM grid storage and shipping puck systems. **Each system will vary as to the type and number of items included with the system.** Potential components include:

- › Cryo-EM Puck
- › Cryo-EM Grid Box Tweezers
- › Cryo-EM Puck Dewar and Insert
- › Cryo-EM Puck Grasping Tongs
- › Bent Cryo Tongs For Puck Manipulation
- › Screw Driver for Puck Lid Lock
- › Cryo-EM Shipping Cane
- › Cryo-EM Storage Cane
- › Barcode Scanner
- › Choice of Puck Color
- › Cryo-EM Puck and Storage Cane Log Sheets
- › Dryshipper with Case Storage Dewar with Rolling Base