

**SPINE Puck Transfer Tool**

for transfer of SPINE sample holders

Cat. No.	Amount
CC-CSM003-0005A	1 pc.

**Description:**

This tool is designed to help loading and unloading SPINE pucks with vials and sample holders. It is intended to sit in a liquid nitrogen bowl such as foam dewars and is designed to elevate the vials/sample holders from their position in the puck for easy removal and transfer to or from another holder. Without this tool, the magnetic ring on the vial keeps the samples stuck inside the puck, rendering the unloading operation very difficult to achieve.

For general laboratory use.**Shipping:** shipped at ambient temperature**Storage Conditions:** store at ambient temperature**Shelf Life:** n/a