

40 µM CELL STRAINER

Catalog # 27305

50 strainers/case

PRODUCT DESCRIPTION:

Sterile 40 µm nylon cell strainer. Useful for obtaining single cell suspensions.

STABILITY / STORAGE:

Store at room temperature.

DIRECTIONS FOR USE:

Unwrap cell strainer unit, fit into a 50-mL centrifuge tube, and filter cells through the strainer.
For single-use only.

THIS PRODUCT IS FOR RESEARCH USE ONLY.