

Catalog #: AS-10134

AnaSpec, Inc.

Product Information Sheet

Seals/13mm, Opentop



For sealing protected amino acid cartridges used in ABI synthesizers.