

ReadiPrep™ Streptavidin Magnetic Beads *3 µm, 1.0%*

Catalog Number: V105080, V105081

Unit Size: 1 mL, 5 mL

Product Details

Storage Conditions	Refrigerated (2-8 °C)
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	N/A

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Streptavidin magnetic beads capture biotinylated biomolecules, enabling separation of antibodies, proteins, antigens, nucleic acids, cells and more.