

**ReadiPrep™ Streptavidin Magnetic Beads \*1 µm, 1.0%\***

Catalog Number: V105070, V105071

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.