

**ReadiPrep™ Streptavidin Magnetic Beads \*0.6 µm, 1.0%\***

Catalog Number: V105060, V105061

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.