

Tosyl-Activated Magnetic Beads *3 µm, 2.5%*

Catalog Number: V105090

Unit Size: 2 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

Tosyl-activated magnetic beads are useful for protein isolation due to the ability of surface groups to covalently couple with amino or sulfhydryl moieties on target biomolecules, such as proteins, antibodies, and glycoproteins. This allows for rapid immunoprecipitation (IP) of proteins and protein complexes.

These magnetic beads are supplied as monodisperse particles (3 µm in diameter) in a 2.5% (25 mg / mL) solution.