

Latex Microsphere *150 nm, White, COOH, 10%*

Catalog Number: V105030

Unit Size: 10 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

Latex microspheres, or latex beads, are colloidal particles composed of amorphous polymer.

Our latex beads are offered in sizes ranging from 80 nm to 300 nm. They are manufactured using high-quality polystyrene and surface coated with carboxyl groups. Additionally, stable, scalable production methods ensure high uniformity in particle size and minimizes batch-to-batch variation, enabling reproducibility in downstream development.

The presence of surface carboxyl groups allows bioconjugation with amine groups on biomolecules- such as peptides, proteins and nucleic acids- through EDC chemistry. This is useful for the development of research tests, for example, solid phase immunoassays.

Density	1.05 g/cm ³
Dispersion medium	Deionized water
Refractive index	1.59 (589 nm, 25 °C)
Functional group	COOH
Bead concentration	10% (100 mg / mL)