



Certificate of Analysis for M-PVA SAV1 Magnetic Beads

Product name	M-PVA SAV1 Magnetic Beads
Cat. no.	CMG-227
Lot no.	SAV1-
Manufacturing date	yyyy-mm-dd
Suspension medium	PBS-Buffer (pH 7.0), 0.05% NaN ₃
Functionalization	Streptavidin
Storage	At +2 to +8 °C
Stability	yyyy-mm-dd (2 years from production date)

M-PVA SAV1 Magnetic Beads

Parameter	Acceptance Criteria	Result
Particle concentration	25.0 ± 1.0 mg/ml	Accord. specification
Particle size	> 90.0% of all particles in the range of 0.60 - 1.25 µm	Accord. specification
Binding properties	> 600 pmol biotinylated OdT ₃₀ per g Beads	Accord. specification
Magnetite content	50 - 60% of total dry weight	Accord. specification
Sterility	No growth of microorganisms visible after 7 days on agar plate	Accord. specification

This Certificate of Analysis has been produced electronically and is valid without a signature. For questions regarding this Certificate of Analysis please contact support.chemagen@revvity.com.