



## Certificate of Analysis for M-PVA OdT2 Magnetic Beads

Product name	<b>M-PVA OdT2 Magnetic Beads</b>
Cat. no.	CMG-231
Lot no.	OdT2-01xxxxx
Manufacturing date	yyyy-mm-dd
Suspension medium	Water
Functionalization	Oligo(dT) <sub>30</sub>
Storage	At +2 to +8 °C
Stability	yyyy-mm-dd (2 years from production date)

### M-PVA OdT2 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 0.80 - 3.0 µm	Accord. specification
Binding properties	> 45 nmol OdA <sub>30</sub> per g Beads	Accord. specification
Magnetite content	50 - 60% of total dry weight	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](mailto:support.chemagen@revvity.com).**