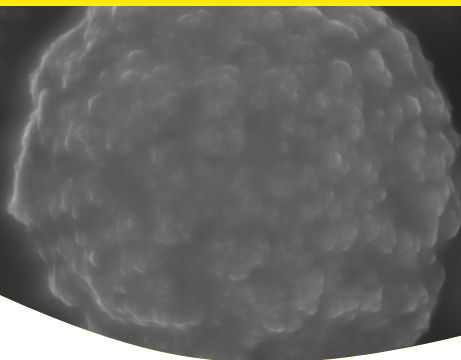




chemagic™ M-PVA Magnetic Beads, where innovation begins.

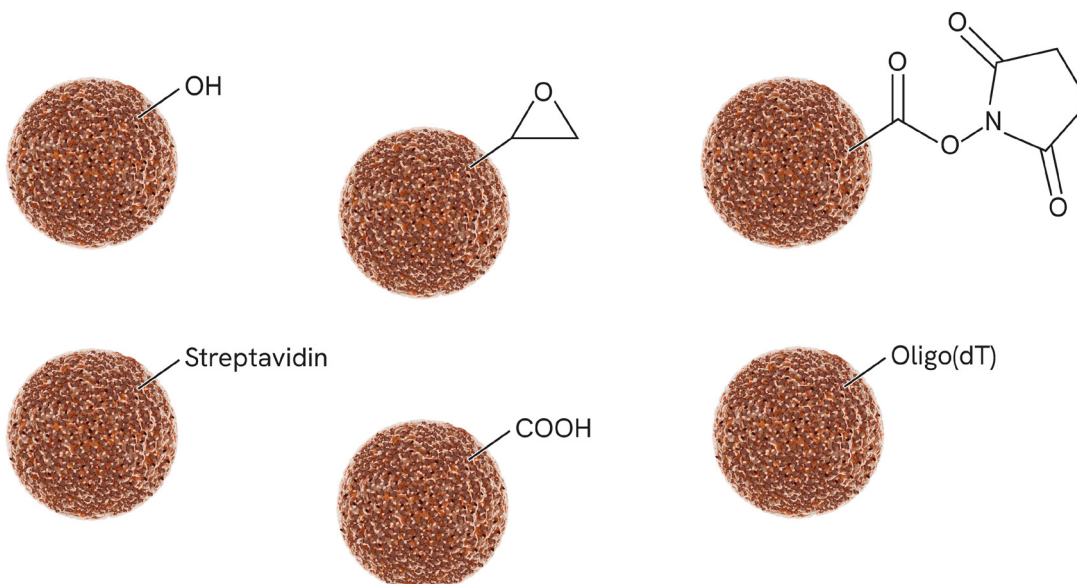


One bead platform, endless possibilities.

From unlimited options for your basic research, over tailored beads for routine assays to ready-to-use applications to complete your product portfolio.

Functionalized magnetic beads

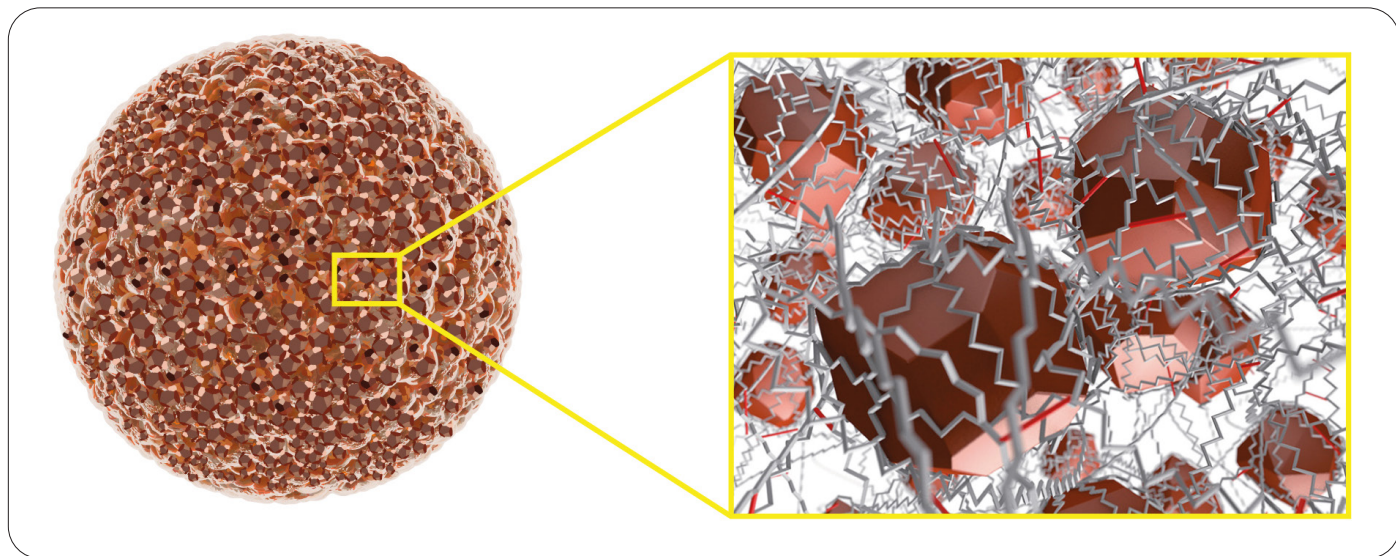
- **Hydroxylation:** Versatile starting point for a wide range of further surface modifications
- **Carboxylation:** Default option for high quality nucleic acid purification, ideal for activation and coupling with EDC/NHS chemistry
- **Epoxidation:** Direct coupling of biomolecules with free amine and thiol groups, supporting your proteome analysis
- **NHS-activation:** Ready-to-couple with proteins and other amine containing ligands, efficient, one-step ligand conjugation
- **Streptavidin:** Target capture of biotinylated probes, high affinity CLIA assay option
- **Oligo(dT):** Specific mRNA capture and enrichment or depletion



Magnetic beads for research & development

Versatile and flexible: discover what's possible

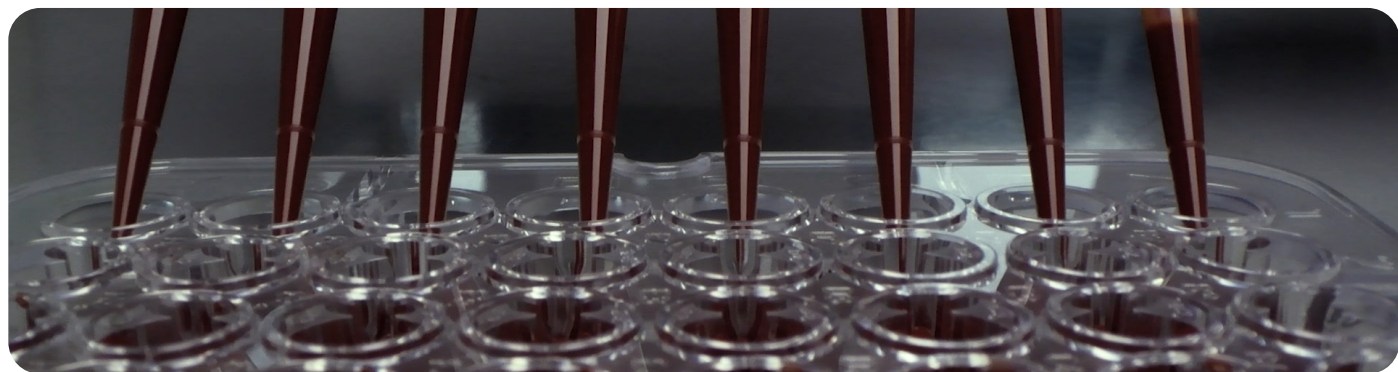
- Cross-linked polyvinyl alcohol matrix entrapping magnetite: Stable and rigid structure with hydrophilic surface
- Superparamagnetic: No residual magnetism, beads redisperse easily after separation
- High magnetite content over 50%
- Two size ranges: 0.6–1.25 μm & 0.8–3.0 μm with tunable binding capacities
- Robustness: Highly tolerable to process solutions
- Stable at +2°C to + 25°C (exceptions apply)



| Cross-linked polyvinyl alcohol matrix entrapping magnetite, stable and rigid structure with hydrophilic surface.

Versatile and flexible: discover what's possible

- Plug-and-play beads for routine diagnostics and research
- Superparamagnetism for efficient working
- Easy pull-down in all sample matrices
- Two bead size ranges (0.6–1.25 μm and 0.8–3.0 μm) with high active surface
- Reliable chemical and mechanical stability
- High tolerance to temperature shifts*, autoclavable



Magnetic bead solutions for biotech, pharma, and life science

Seamless integration into your workflow

- Instrument friendly: Short run times in your instrument
- Compatible with large volume samples
- Application dedicated bead size
- Physical strengths for automation
- Storable in water*, easy shipment, colerant to lyophilization
- Stringent QC release criteria

**not applicable for all magnetic beads*



Trusted expertise. Proven quality. Dedicated partnership.

With **over 25 years of experience** in bead production and functionalization, we bring deep expertise and a proven track record to every partnership.

Our commitment to quality is backed by both **ISO 9001** and **ISO 13485** certifications, providing that our product meets the standards your applications demand.

As a **single-source supplier**, we simplify your supply chain by providing what you need under one roof — from research and application support to full-scale production. With our established production capabilities, we are equipped to scale with your needs.

Our **flexible kit configurations** allow you to tailor solutions to your specific requirements, while our **1:1 dedicated OEM support** provides you have a committed partner by your side — from development through delivery.



Part number	M-PVA code	Description	Particle size (µm)	Concentration (mg/mL)	Unit (mg)	Storage (°C)	Shelf life (months)
Hydroxylation: versatile tool for a wide range of further surface modification							
CMG-201	M-PVA 012	Hydroxylated M-PVA Magnetic Beads	0.8 – 3.0	50	500	+2 to +25	36
Carboxylation: default option for high quality nucleic acid purification							
CMG-203	M-PVA C11	Medium carboxylated M-PVA Magnetic Beads	0.6 – 1.25	50	500	+2 to +25	36
CMG-206	M-PVA C21	Highly carboxylated M-PVA Magnetic Beads	0.6 – 1.25	50	500	+2 to +25	36
CMG-207	M-PVA C22	Highly carboxylated M-PVA Magnetic Beads	0.8 – 3.0	50	500	+2 to +25	36
Epoxidation: direct coupling of biomolecules with free amine and thiol groups							
CMG-216	M-PVA E02	Epoxydated M-PVA Magnetic Beads	0.8 – 3.0	50	500	+2 to +8	6
NHS-activation: ready-to-couple with proteins and other amine containing ligands							
CMG-221	M-PVA Ak11	NHS activated M-PVA Magnetic Beads	0.6 – 1.25	25	50	+2 to +8	3
CMG-222	M-PVA Ak12	NHS activated M-PVA Magnetic Beads	0.8 – 3.0	25	50	+2 to +8	3
CMG-237	M-PVA Ak11	NHS activated M-PVA Magnetic Beads	0.6 – 1.25	25	500	+2 to +8	3
CMG-238	M-PVA Ak12	NHS activated M-PVA Magnetic Beads	0.8 – 3.0	25	500	+2 to +8	3
Streptavidin: target capture of biotinylated probes							
CMG-227	M-PVA SAV1	Streptavidin coated M-PVA Magnetic Beads	0.6 – 1.25	25	50	+2 to +8	24
CMG-228	M-PVA SAV2	Streptavidin coated M-PVA Magnetic Beads	0.8 – 3.0	25	50	+2 to +8	24
Oligo(dT): specific binding of mRNA							
CMG-231	M-PVA OdT2	Oligo(dT) coated M-PVA Magnetic Beads	0.8 – 3.0	25	50	+2 to +8	24

Need something beyond the catalog?

Your research is unique, your beads can be too.

If your project requires specific particle properties, a custom ligand, or a conjugation chemistry not listed here, our team of experts is ready to collaborate. We support you from concept to delivery.



Learn more!

**Scan or click
for more information**

revvity