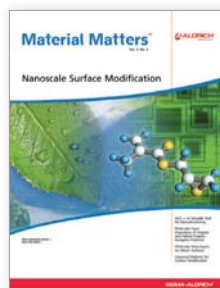


Innovative Materials from Sigma-Aldrich

You may be familiar with Sigma-Aldrich research chemicals for synthesis, but did you know we also offer innovative materials for your research needs into newly emerging technologies?

For a free copy of our *Materials Science Catalogue* please visit sigma-aldrich.com/mscatalogm



...and whilst you are online why not subscribe to complimentary copies of *Material Matters*, a quarterly publication dedicated to addressing your research needs, featuring technical reviews, application notes and relevant services.

sigma-aldrich.com/mmca

For a quotation on any of our products please contact us at uksalesupport@sial.com or for Technical Service advice email us via eurtechserv@sial.com.

Alternative Energy — For your research into clean sources of energy, including materials for hydrogen storage

Metal and Ceramic Science — Offering high purity metals, metal oxides and metal salts as well as metal alloys, ceramics and binary and ternary metal compounds

Micro/NanoElectronics — Products designed for application in chemical vapour deposition (CVD), chemical solution deposition (Sol-Gel), atomic layer deposition (ALD), lithography and nanopatterning, molecular self-assembly and many others. Electronic chemicals and electronic materials are also available.

Nanomaterials — Addressing your nanotechnology needs from research into drug delivery through to nanodevices. We have inorganic nanopowders, nanoparticle dispersions and surface functionalised nanoparticles. Carbon nanotubes, fullerenes, dendrimers, dendrons, hyperbranched polymers, nanoclays and quantum dot nanocrystals are included too.

Organic Electronics and Photonics — Consisting of three groups, organic electronics, photonics and liquid crystals

Polymer Science — Spans monomers, polymers including biodegradable polymers, polymer additives and polymerization initiators.

Products for research into the above areas can be viewed in our Materials Science Catalogue and as our range is continually expanding, many more are listed on our website sigma-aldrich.com/matsci.

Sigma-Aldrich Company Ltd.
The Old Brickyard
New Road
Gillingham
Dorset SP8 4XT

Order/Customer Service Phone 0800 71 71 81 • Fax 0800 378785
ukorders@sial.com

Technical Service 0800 272572 • sigma-aldrich.com/techservice

Development/Bulk Manufacturing Inquiries SAFC® 01202 712305

*Accelerating Customers' Success
through Innovation and
Leadership in Life Science,
High Technology and Service*