

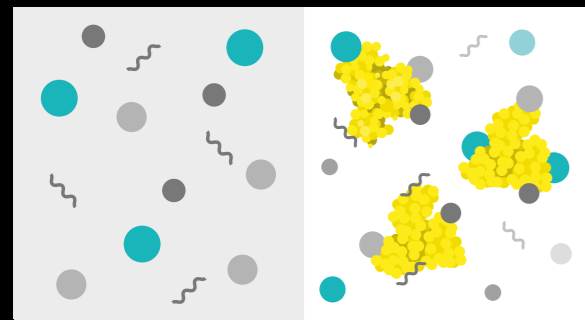
Recombunin[®] Recombinant Human Albumin

Your Multifunctional Tool From Concept To Manufacturing and Distribution

Recombunin[®] recombinant human albumin (rHA) is a high-quality animal- and human-origin-free rHA. It meets cGMP manufacturing standards and has USP-NF status, and its stabilizing properties have been validated through years of use in multiple life science products.

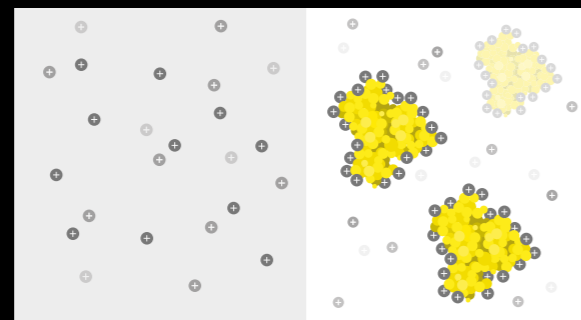
Mechanisms of Action | How Recombunin[®] Delivers Value to Your Process

Biological Mechanisms of Action



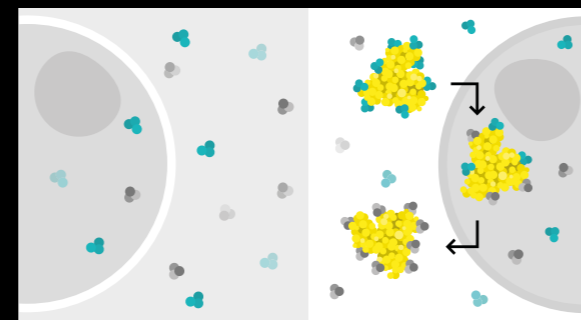
Nutrient Carrier

Binds and transports nutrients but is itself a rich nutrient source for cells



pH Buffer

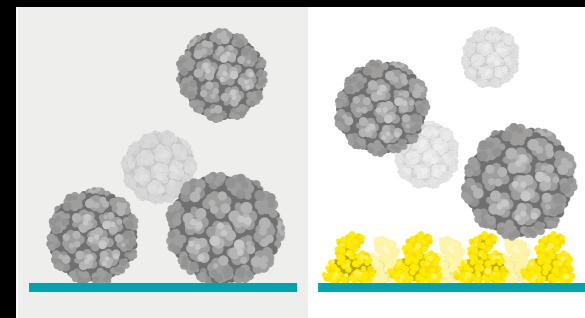
Binds to hydrogen ions and buffers pH



Transportation

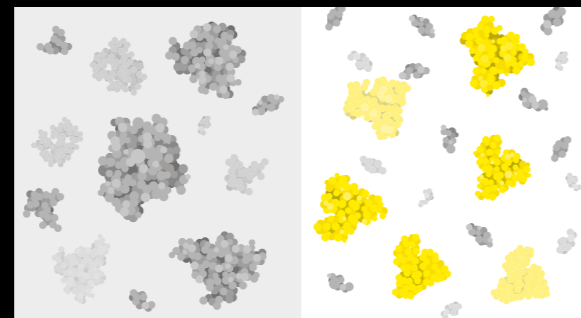
Forms complexes with metals and molecular entities transporting these into and out of cells

Physicochemical Mechanisms of Action



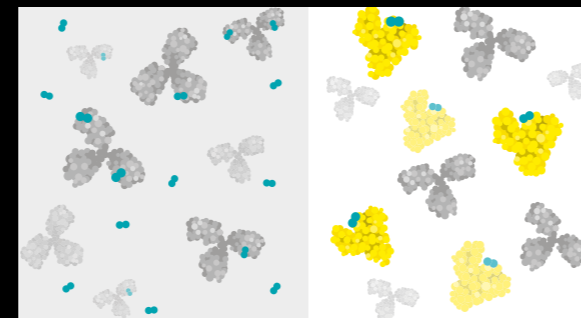
Adsorption

Forms a single monolayer to prevent adsorption



Aggregation

Prevents aggregation and disperses uniformly in a solution



Oxidation

Protects formulations from oxidative stress

Supported Applications

Recombunin[®] can confer a range of benefits when used in different applications across several modalities; supporting products from the laboratory bench to clinical application.



Stem Cell Therapies



Gene Modified Cell Therapies



Viral Therapies and Vaccines



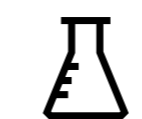
Protein Based Therapies



Gene Therapies



Medical Devices



Diagnostics and Reagents

Designed by Nature, Enhanced by Sartorius

Quality

Manufactured to cGMP standards. Consistent, pure, and animal- and human-origin-free.

Performance

Offers high performance backed up by years of use in preclinical, clinical, and marketed biopharmaceuticals and medical devices.

Regulatory

Comes with comprehensive regulatory documentation and support that leverages our near four decades of expertise.

SARTORIUS