Recombumin[®] Recombinant Human Albumin | Medical Device Applications

The physicochemical properties of Recombumin[®] recombinant human albumin (rHA) offer a solution to biocompatibility challenges in medical device manufacturing, such as adsorption and biological-mechanical interfaces.

For developers working on the next generation of advanced medical devices, the biocompatibility of coating is crucial. Albumin's natural properties and ubiquitous presence in the body make Recombumin[®] rHA the ideal choice for developers.

Hydrophobic and Ionic **Properties Allow** Recombumin[®] to Coat a Range of Surfaces

- Stops surface adsorption to optimize yield while controlling dose reliability and precision
- Forms a single monolayer that can coat 1m² of surface area using just 1-2 mg of rHA
- Offers a biocompatible human- and animalorigin-free surface coating



Creating an Anti-adsorptive and Biocompatible Coating for ECMO Blood Oxygenation Devices.

Chalice Medical Ltd. is a UK-based medical device manufacturer making high-quality tubing packs and oxygenators in ISO-7clean room conditions.







Case Study



Their extracorporeal membrane oxygenation (ECMO) devices support direct oxygenation of human blood.

Chalice needed a biocompatible, antiadsorptive surface coating.

Recombumin[®] rHA was selected to be a

component of their surface coating for ECMO devices.



Your Benefits

- Chemically inert
- USP-NF status
- cGMP compliant
- Comprehensive regulatory documentation
- Biologically stable
- Solvent-free
- Nontoxic
- Biocompatible

SVIDENS