# SVISCISAS

### **Success Story**

## Enabling Rapid Vaccine Production Scale-up

#### **Case Profile**

Company Type: Small | Startup biotech, CDMO or biopharma

Related Molecule: Vaccines

Process Steps: Drug Substance Storage, Shipping, and Supply Chains

A small biopharma company needed a quick solution on shipping substances to various locations and the potential to scale up to larger volumes going through clinical phases and bulk production in a very short timeframe.

#### **Customer Challenge**

- Rapid development and commercialization of vaccine
- Transportation of drug substances between sites
- Upscaling of production required additional personnel that needed to be trained in a short timeframe

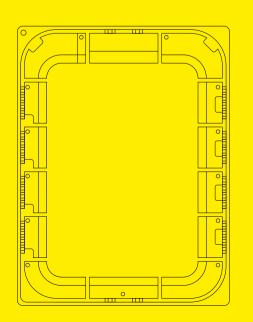
#### **Provided Solution**

- Sartorius provided solution to ship frozen drug substances safely and reliably (products in focus: Celsius® FFT with Safecore technology and Celsius® FFT bulk shipper)
- Sartorius enabled quick ramp up of production scales of vaccine and provided solutions in close alignment on scale-up needs of customer (immediate response to quantities needed)
- Implementation and operator adoption was supported by the products' ergonomic and easy-to-use design. Additionally, employee training programs were initiated to simplify effective product handling

>100 employees trained in on Celsius<sup>®</sup> FFT within 6 months

3,000% increase of vaccine production within 6 months

50 million vaccine doses produced in 6 months



"Time was a critical factor. Sartorius enabled efficient production scale-up for our customer by responding quickly to their needs, providing expertise and support, and ensuring products were available within a very short timeframe."

**Elisabeth Pfuetzner-Riehn** Sales Representative - General Sales

Contact a Sartorius Specialist or Visit www.sartorius.com/freeze-thaw