

May 14, 2021

Dear Valued Customer:

We are pleased to announce the global expansion of Thermo Fisher Scientific Production Chemicals and Services in Europe through the distribution agreement with Avantor, effective January 1, 2021. The catalog offering of cGMP production chemicals includes specific products from Avantor, one of the world's leading chemical suppliers for biomanufacturing applications. Additional catalog chemical suppliers will be added in the upcoming months as we continue to expand our global footprint. The list below itemizes the specific chemical categories available to our customers through the Avantor distribution agreement.

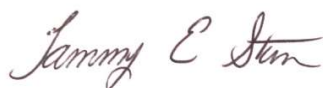
Available Chemical Categories

- Acid/base solutions
- Amino acids
- Biological buffers
- Biological reagents
- Carbohydrates
- Cleaning agents and disinfectants
- Denaturants
- Emulsifiers
- Media supplements
- Minerals
- Salts
- Solvents
- Vitamins

Access to these new catalog chemical products complements our current offering of Process Liquid Preparation Services. The catalog chemical product offering allows the convenience of "off the shelf" products to be purchased, while the Process Liquid Preparation Services is designed to simplify the process liquid and buffer preparation workflow to enable bioprocessing customers to produce life-changing therapeutics and vaccines, faster. Through close collaboration, we work with customers to create fit-for-purpose solutions ranging from ready-to-hydrate dry powder chemicals, ready-to-use process liquids, or a combination of both, including flexible options for chemical grade, format, and container type.

We look forward to serving you with catalog cGMP Production Chemicals and Process Liquid Preparation Services to simplify your bioproduction workflows.

Sincerely,



Tammy Starr
Vice President/General Manager
Thermo Scientific Production Chemicals and Services