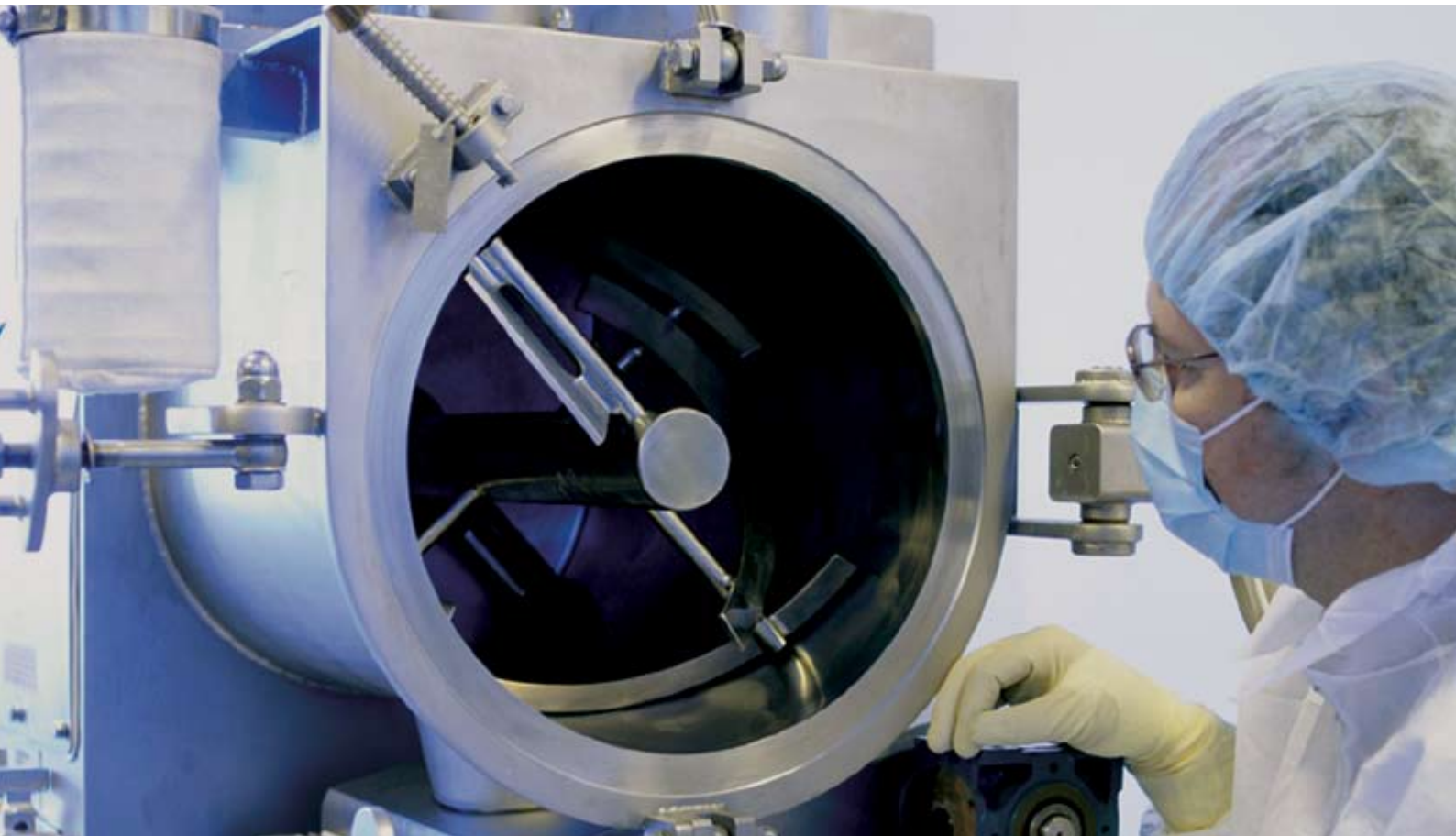


## Powdered Media State-of-the-art Facility Now Open



We are pleased to announce the commissioning of our new powdered media production facility in Walkersville, Maryland. This state-of-the-art, non-animal origin, cGMP manufacturing suite is designed for the efficient production of media batches from 7.25 – 7,290 kg using micronizing mills and ribbon blenders.

The newly renovated suite offers approximately 3,000 square feet of manufacturing space that is temperature and humidity controlled, with continual environmental monitoring. The equipment, designed for continuous operation, includes 2 mills and 4 ribbon blenders, all stainless steel construction, and is equipped for clean in place (CIP) for ease of use.

From pilot to commercial scale, we can produce a product to meet your needs. We invite you to tour or audit our new facility today.



## Batch Sizes for Minimum and Maximum Loads for Blenders

Blender	Minimum Load		Maximum Load	
	kg	liter	kg	liter
50 L	7.25	725	40.5	4,050
300 L	43.5	4,350	243	24,300
2,000 L	290	29,000	1,620	162,000
9,000 L	1,305	130,500	7,290	729,000

Our new powdered media production area is set up with two independent systems:

- The small train includes our small-scale micronizer and 50 L and 300 L blenders. This is a semi-manual system designed to meet the needs of small-scale productions from approximately 725 L to 24,300 L of powdered media. It is manually dispensed and cleaned using a fully validated clean in place (CIP) skid.
- The large train system includes a large-scale micronizer and 2,000 L and 9,000 L ribbon blenders. Bar coded raw materials are sequentially added to the mill and conveyed pneumatically through a completely closed system until final dispensing, preventing contamination of the product and the environment. This train is equipped with spray balls for complete CIP using the validated cleaning skid.

Both systems deliver uniform particle size where  $\geq 95\%$  of particles are  $< 212$  microns, the optimal size for dissolution of dry powdered media formulation.

Lonza powdered media manufacturing offers the following advantages:

- Raw materials of non-animal origin create a regulatory friendly facility
- Consistent particle size distribution results in excellent media solubility
- High-speed impact milling avoids chemical degradation
- Automated CIP technology avoids contamination between batches



## Contact Information

### North America

Customer Service: 800-638-8174  
Technical Service: 800-521-0390  
E-mail: [biotechserv@lonza.com](mailto:biotechserv@lonza.com)  
Online Ordering: [www.lonza.com](http://www.lonza.com)

### Europe

Customer Service: 32 (0) 87 321 611  
Technical Service: 32 (0) 87 321 651  
(German Speaking)  
32 (0) 87 321 626  
(French Speaking)  
E-mail: [techsup.europe@lonza.com](mailto:techsup.europe@lonza.com)  
or [techsup.uk@lonza.com](mailto:techsup.uk@lonza.com)  
Online Ordering: [www.lonza.com](http://www.lonza.com)

### International

Contact your local Lonza Distributor  
Customer Service: 301-898-7025, ext. 2322  
Fax: 301-845-8291  
E-mail: [biotechserv@lonza.com](mailto:biotechserv@lonza.com)

### International Offices

Australia 61 3 9550 0883  
Austria 0800 201 538  
Belgium 32 (0) 87 321 611  
Denmark 45 43 56 74 07  
France 0800 91 19 81  
Germany 0800 182 52 87  
India 91 22 6697 2883  
Italy 39 0363 35 14 70  
The Netherlands 0800 022 4525  
Spain 34 902 531 366  
Switzerland 0800 83 86 20  
United Kingdom 44 118 979 5234

Lonza Walkersville, Inc.  
Walkersville, MD 21793

For further manufacturing and laboratory use. Not for use in diagnostic procedures.

Unless otherwise noted, all trademarks herein are marks of the Lonza Group or its affiliates.

© Copyright 2007, Lonza Walkersville, Inc. All rights reserved.  
FL-POWDER-1 08/07

**We cordially invite you to tour our new powdered media manufacturing facility.**

We are located in Walkersville, MD, convenient to Baltimore, Washington-Reagan National and Dulles airports. Contact us today to arrange a site visit.

Lonza Walkersville, Inc. 800-521-0390

**Lonza**