



Notice of Upcoming Changes to Product Codes and Disease State Product Labels

To better serve you, we are improving our product management system. The following changes will make it easier for you to place orders with us:

1. Code updates for all products, except ADME and ADME service offerings
2. Product label updates for disease state biofluids
3. Standard volume changes to disease state biofluids

Keep reading for additional details about these changes.

Product Codes

Product codes are being standardized and encoded to capture key product defining elements, such as:

- Format
- Diagnosis
- Cell number per vial
- Shipping media
- Charcoal stripping
- Filtration

Note that going forward, codes will be reusable for products with the same alignment.

Disease State Product Labels

Product labels will no longer designate a single disease indication. Instead, they will contain a 2D barcode and the following information:

- Part number
- Lot number
- Sublot
- Volume
- Storage temperature

Additionally, orders will come with a digital file containing a list of specimens with clinically relevant information, such as diagnosis, medications, test results and social demographic information.

Disease State Standard Volume

We are pleased to announce the standardization of all blood-derived disease state biofluids to 1mL aliquots.

This new process offers several advantages, including:

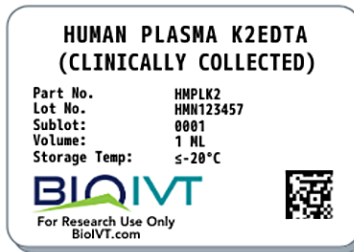
- Increased standardized protocols for collection and processing
- The ability to view product inventory online with pricing and purchasing options

All of these adjustments are ongoing, and you will begin to see changes implemented in January. We realize that you may require additional support to update your systems. As always, our client service representatives are available to assist you. Please do not hesitate to reach out to us at customerservice@bioivt.com or cseurope@bioivt.com.

Product Code Transition Examples

Product Description	Current Product Code	Ex. New Product Code
Human Plasma K2EDTA (normal)	HUMANPLK2UZN	HUMANPLK2-#####
Human Plasma K2EDTA Lung Cancer	HMPLK2LNGCA or HMPL-LNGCA2	HUMANPLK2-#####
Human LEUKOMAX® Mobilized Leukopak 5-day Neupogen/1 day Mozobil, collected on day 5	HUMANLMX5NM1	HUMANMLMX515-#####
Beagle Whole Blood Ultrafiltered, Charcoal Stripped 3x	CAN00WBU2103	CAN00WB-#####
CD-1 (ICR) Mouse PBMC cells, whole blood, 5x10 ⁶ cells/vial, mixed gender, pooled from 10 mice	MSE00-PBMC-P-201699	MSE00PBMC-#####
Human Lung Cancer Adenocarcinoma FFPE	HMLNG-ADENO1	HUMANLNG-#####

New Product Label Example



Follow us on social media:



BioIVT

North America & Asia Pacific
PO Box 770, Hicksville, NY, 11802-0770
T 516-483-1196
customerservice@bioivt.com

Europe, Middle East & Africa
T + 44 (0) 1444 707333
cseurope@bioivt.com

You received this email because you are subscribed to Marketing Information from BioIVT.

Update your [email preferences](#) to choose the types of emails you receive.

[Unsubscribe from all future emails](#)

[Contact Us](#) | [Privacy Policy](#) | [Terms & Conditions](#)