

Product Change Announcement

Echelon has changed the PIP₃ Mass Strip kit (K-2400). Changes include:

1. A new format of PIP₃ Mass Strip with PI and 6 other PIP_n control spots, 8 PIP₃ standard spots, and room for users to spot 8 unknown samples.
2. A new PIP₃ detector protein in lyophilized form. Three vials are supplied for three separate assays.
3. Different secondary and tertiary detectors.
4. Revised technical data sheet and protocol.

The new PIP₃ detector protein has proved to be more sensitive and specific for PIP₃ when compared to the original PIP₃ detector. **The product itself is identical in all other aspects of performance and specifications.** Customers should verify that they are following the correct technical data sheet/protocol before using the product. Echelon regrets any confusion this change might cause. Our intention is to provide the highest quality research products to our customers. From time to time as we continue to improve our products, Echelon will use “Product Change Announcements” to communicate these changes.

These changes will affect kits produced after 11/27/2007 (with lot numbers greater than XX112707). For technical questions please contact Kim Bruner at kbruner@echelon-inc.com or call 801-588-0455 ext. 305.