

Product Change Announcement

Echelon has changed the Class III PI3-K ELISA kit (K-3000). Changes include:

1. PI(3)P Detector (K-3005) has been replaced by lyophilized form of the same protein with cat# K-3305. K-3305 has proved to be more stable than K-3005. Four vials of K-3305 are supplied for up to 4 separate assays.
2. Secondary Detector (K-3006) and Tertiary Detector (K-3007) have been replaced by a single HRP Conjugate (K-2306). This change has simplified the assay and increased the signal window.
3. Following reagents have been added to the kit to make it more user friendly:
 - 10x TBS-T buffer (K-TBST)
 - 100 mM EDTA (K-EDTA)
 - 1 N H₂SO₄ Stop Solution (K-STOPt)
4. Revised technical data sheet and protocol.

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. 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 01/01/2008 (with lot numbers greater than XXX010108). For technical questions please contact Xin Morrow at xchen@echelon-inc.com or call 801-588-0455 ext. 375.