



P.O. Box 58537
Salt Lake City, Ut 84108
Telephone 801-588-0455
Fax 801-588-0497
Email: echelon@echelon-inc.com
http://www.echelon-inc.com

Certificate of Analysis

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Product Name: PI 3-K Fluorescence Polarization Activity Assay

Product Number: K-1100

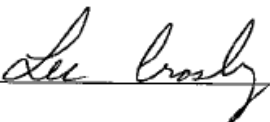
Lot Number: AB102009

Product Description: FP kit including reaction plate, PI(4,5)P₂ Substrate, PI(3,4,5)P₃ Standard, PI(3,4,5)P₃ detector, Red Fluorescent Probe, PI(3,4,5)P₃ Detector Diluent, Probe Diluent, and black, non-binding 384 well plate.

Quality Assay: Fluorescence Polarization

Quality Parameters	Required	Actual	Passes
A. R ²	>0.95	0.97	✓
B. Z'	>0.4	0.85	✓
C. Specificity	>2	2.9	✓
D. Sensitivity	1±0.2	1.18	✓
E. Window	>2	2.5	✓

Passes All Functional Specifications



Date 10/21/2009

Lee Crosby
Quality Manager