

# Acyl Labeled PIP's



**Echelon's  
back in the  
game!**

**-Available August 1, 2008-  
Echelon offers labeled  
analogs for all 8 PIP head  
groups with biotin and  
fluorescent tags.**



**echelon**  
www.echelon-inc.com