



Echelon Biosciences Inc.
675 Arapeen Dr., Ste. 302
Salt Lake City, UT 84108
Phone: 801.588.0455
Fax: 801.588.0497

ACYL LABELED PIPS

Available August 1, 2008

Headgroup	Cat #	Biotin	Label	
			BODIPY [®] FL	BODIPY [®] TMR
Phosphatidylinositol (PtdIns)	Cat #	C-00B6	C-00F6	C-00M6
Phosphatidylinositol 3-phosphate (PtdIns(3)P)	Cat #	C-03B6	C-03F6	C-03M6
Phosphatidylinositol 4-phosphate (PtdIns(4)P)	Cat #	C-04B6	C-04F6	C-04M6
Phosphatidylinositol 5-phosphate (PtdIns(5)P)	Cat #	C-05B6	C-05F6	C-05M6
Phosphatidylinositol 3,4-bisphosphate (PtdIns(3,4)P ₂)	Cat #	C-34B6	C-34F6	C-34M6
Phosphatidylinositol 3,5-bisphosphate (PtdIns(3,5)P ₂)	Cat #	C-35B6	C-35F6	C-35M6
Phosphatidylinositol 4,5-bisphosphate (PtdIns(4,5)P ₂)	Cat #	C-45B6	C-45F6	C-45M6
Phosphatidylinositol 3,4,5-trisphosphate (PtdIns(3,4,5)P ₃)	Cat #	C-39B6	C-39F6	C-39M6