



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

New Product Update:

June 16, 2008

Assay Services (send us your sample and we'll run the assay):

- Available 7-1-08 -

Echelon catalog #	Product Name
T-1900	Sphingosine 1 Phosphate ELISA
T-2800	Lysophosphatidic Acid ELISA
T-1200	Hyaluronic Acid ELISA
T-1700	Collagen IV ELISA
T-31SM	Sphingomyelin Assay
T-31PC	Phosphatidylcholine Assay
T-1800	Sphingomyelinase (neutral) Assay
T-3200	Sphingomyelinase (acidic) Assay
T-2300	PTEN Assay
T-1400	SHIP2 Assay
T-0LBT	Lipid Binding Testing Service

Echelon catalog #	Product Name
M-1000	PIP Microarray Plate
K-3200	Acidic Sphingomyelinase Assay
Z-R015	Anti-HVps34 Antibody for immunoprecipitation
Z-R016	Anti-HVps34 Antibody for western blot
G-0302	PI(3)P Grip
Z-DGK α	Anti-diacylglycerol kinase α polyclonal antibodies
Z-DGK ζ	Anti-diacylglycerol kinase ζ polyclonal antibodies
S-1000	Sphingosine
S-2000	Sphingosine 1 Phosphate
S-200B	Sphingosine 1 Phosphate Biotin
C-39F16a	Bodipy FL PI(3,4,5)P ₃
B-0720	FTY-720
B-0222	Wortmannin
B-0312	OSU-03012 – PDK1 inhibitor
B-0021	Resveratrol
B-0022	Trismethoxyresveratrol
L-7416	BrP-LPA (LysoPLD Autotaxin Inhibitor)
I-M052	4-Diphosphocytidyl-2-C-Methyl-D-erythritol (CDP-ME)
P-Bcer	Ceramide Beads
P-BLBPA	Lysobisphosphatidic Acid Beads
P-B0SM	Sphingomyelin Beads

NEW PRODUCT CATEGORY – HEME PATHWAY INTERMEDIATES

Echelon catalog #	Product Name
F-H010	Hemin
F-H020	Porphobilinogen
F-H030	Uroporphyrin III Dihydrochloride
F-H040	Uroporphyrin I Dihydrochloride
F-H050	Protoporphyrin IX
F-H100	Biliverdin Hydrochloride
F-H110	Biliverdin Dimethylester
F-H120	Bilirubin
F-H130	Bilirubin Conjugate
F-H140	Urobilin Hydrochloride
F-H150	Stercobilin Hydrochloride

Misc. Research Chemicals

Echelon catalog #	Product Name
F-H060	Zn (II) Protoporphyrin IX (Inhibitor of Heme Oxygenase)
F-H070	Cu (II) Protoporphyrin IX (negative control of inhibitor)
F-H080	Sn (IV) Protoporphyrin IX Dichloride (Inhibitor of Heme Oxygenase)
F-H090	Zn (II) Deuteroporphyrin IX 2,4 bis ethyleneglycol (Inhibitor of Heme Oxygenase)
F-0030	Monobromobimane (Thiol Reaction Probe)
F-0040	Dibromobimane (Thiol Reactive Probe)
F-0010	Dimethyloxalylglycine (DMOG) (HIF-1 α stabilizer)
F-0020	N-oxalyl glycine (NOG) (alpha ketoglutarate analogue)