

Cayman Chemical

- Fatty Acyls (Eicosanoids)
- Glycerolipids
- Sphingolipids
- Glycerophospholipids
- Sterol Lipids
- Prenol Lipids

