

## Examples of Agonists that Induce Phosphatidylcholine Breakdown

Vasopressin  
Phorbol esters  
Thrombin  
Acetylcholine  
N-formyl-Met-Leu-Phe  
Ca<sup>2+</sup> ionophores  
EGF  
PDGF  
IL1, IL2, IL3