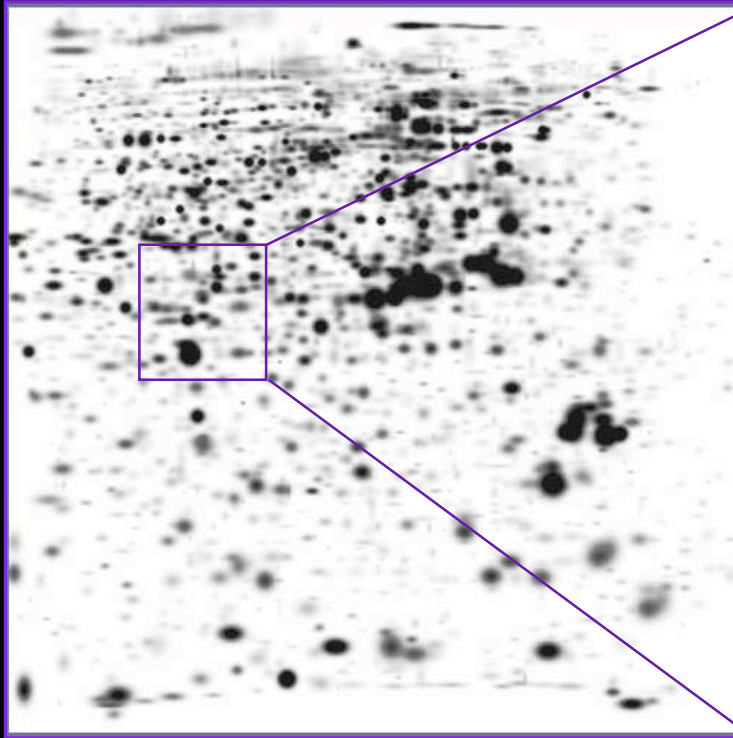
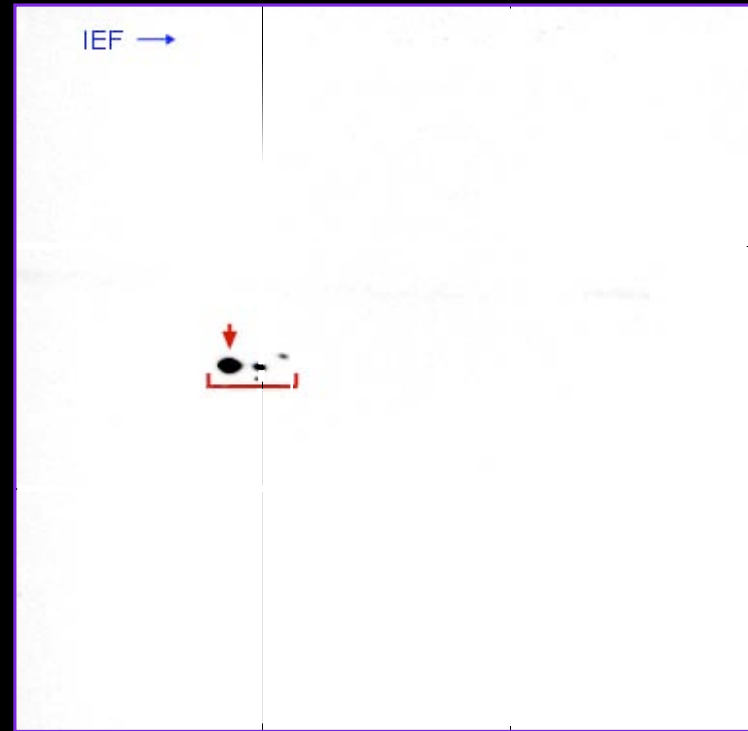


2D Gel Electrophoresis - Immunoblot

- Protein kinases are in low abundance and usually not detectable by silver staining



Silver-stained



Immunoblot with ERK1/2 Ab