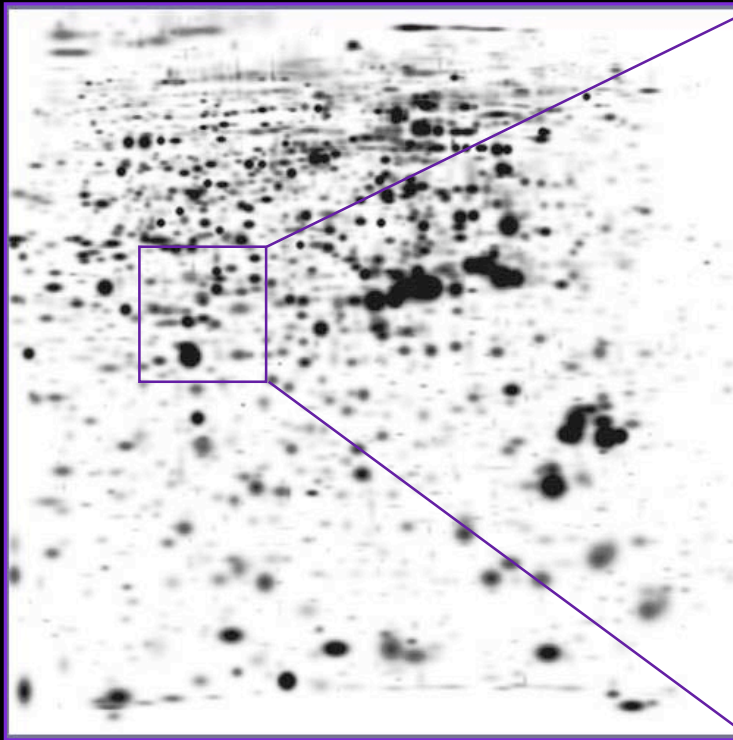


2D Gel Electrophoresis -Silver Stained

Human keratinocyte proteome analysis



Silver-stained

