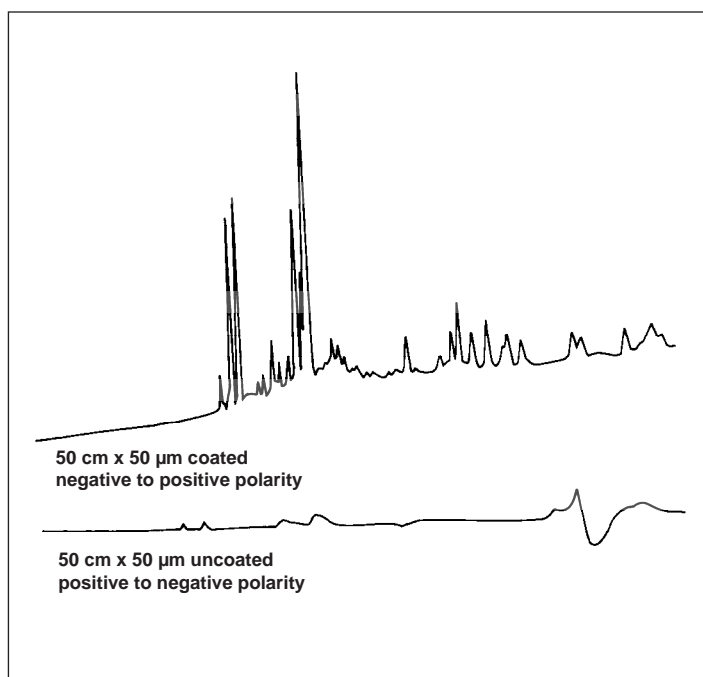


## BSA Tryptic Peptides - A Comparison of Separations in Coated and Uncoated Capillaries

Bio-Rad's patented coatings enhance protein separations at alkaline pH. In addition to reducing solute adsorption, the coating greatly diminishes electroendosmosis. When an uncoated tube is used under the same conditions, most of the peptide components strongly adsorb to the capillary surface. This is illustrated below with capillary electrophoresis of a tryptic digest of bovine serum albumin (BSA) in both coated and uncoated capillaries. Peptides are loaded and separated by electrophoretic mobility on the coated tube, but loaded and separated by electrophoresis and electroendosmosis on the uncoated tube, necessitating reverse polarity.



<b>Capillary:</b>	50 cm x 50 µm, coated and uncoated
<b>Buffer:</b>	0.1 M potassium borate, pH 8.5, linear polymers, zwitterion additives
<b>Load conditions:</b>	8 kV, 8 seconds
<b>Run conditions:</b>	8 kV, constant voltage, negative polarity (coated) and positive polarity (uncoated)
<b>Detection:</b>	UV, 200 nm, 0.1 AUFS