

The outline on this column is made up of multiple vertical segments, two for each text field (left, right). They stretch with the corresponding text field. This solution has the disadvantage that it is impossible to suppress a text field space when the section overflows to a new page. Check the top of the second page to see that.

In order to minimize the space occupied by the text fields even when they do not print, their height was set to zero. There still seems to be some problems with the bottom line that would sometimes print and sometimes not, when the section breaks, depending on the remaining space.

This solution is similar with the previous. The left and right borders are made up of a single vertical line that stretches with the group represented by the three text fields. The page break behavior is almost the same.

This seems to be the best solution for creating an outline for the three text fields. It is made up of a single rectangle that stretches with the group that contains the three text fields. The page breaks seem to work very well.

| Page Header | | | Page 1 |
|--|--|--|--|
| <div>This is a FIRST long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a SECOND long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> | <div>This is a FIRST long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a SECOND long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> | <div>This is a FIRST long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a SECOND long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> | <div>This is a FIRST long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a SECOND long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> |
| <div>This is a FIRST long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a SECOND long chunk of text</div> | <div>This is a FIRST long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a SECOND long chunk of text</div> | <div>This is a FIRST long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a SECOND long chunk of text</div> | <div>This is a FIRST long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.</div> <div>This is a SECOND long chunk of text</div> |
| Page Footer | | | Page 1 |

This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a THIRD
chunk of text that
will cause the text
field to stretch

This is a THIRD
chunk of text that
will cause the text
field to stretch

This is a THIRD
chunk of text that
will cause the text
field to stretch

This is a THIRD
chunk of text that
will cause the text
field to stretch

This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a THIRD chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a SECOND long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a SECOND long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a SECOND long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

This is a SECOND long chunk of text that will cause the text field to stretch outside its defined height and force other elements to move downwards.

