

This is a **static text** element containing styled text. **Styled text** elements are introduced by setting the *isStyledText* boolean attribute available for the **textElement** tag to *true* and by formatting the text content using nested **style** tags and simple HTML tags.

This is a text field element that contains styled text displaying the text-only version of the ^JasperReports logo.

Here is the list with the attributes supported by the **style** tag used to format the text content inside a styled text element:

- fontName
- size
- isBold
- isItalic*
- isUnderline
- ~~isStrikeThrough~~
- pdfFontName
- pdfEncoding
- isPdfEmbedded
- forecolor
- backcolor

Here is the list with the supported HTML tags :

```
<font size="12">  
<font color="red">  
<font face="Comic Sans MS">  
bold  
italic  
underlined  
an empty line follows(<br/>):
```

A bulleted list:

- item 1

ampersand & less < greater > quote "
ampersand & less < greater > quote "

	1	
	2	
	3	
	2	
	1	