

```
<textField><reportElement x="0" y="50" width="200"
height="90"/><textElement><font fontName="Arial" size="14"
pdfFontName="ARIAL.TTF" pdfEncoding="Identity-H"
isPdfEmbedded="true"/></textElement><textFieldExpression>$P{GreekText
} + "    " + $P{CyrillicText}</textFieldExpression></textField>
```

```
<textField><reportElement x="0" y="150" width="200"
height="90"/><textElement><font fontName="Arial" size="14" isBold="true"
pdfFontName="ARIALBD.TTF" pdfEncoding="Identity-H"
isPdfEmbedded="true"/></textElement><textFieldExpression>$P{GreekText
} + "    " + $P{CyrillicText}</textFieldExpression></textField>
```

```
<textField><reportElement x="0" y="250" width="200"
height="90"/><textElement><font fontName="Arial" size="14" isItalic="true"
isUnderline="true" pdfFontName="ARIALI.TTF" pdfEncoding="Identity-H"
isPdfEmbedded="true"/></textElement><textFieldExpression>$P{GreekText
} + "    " + $P{CyrillicText}</textFieldExpression></textField>
```

```
<textField><reportElement x="0" y="350" width="200"
height="90"/><textElement><font fontName="Arial" size="14" isBold="true"
isItalic="true" isUnderline="true" pdfFontName="ARIALBI.TTF"
pdfEncoding="Identity-H"
isPdfEmbedded="true"/></textElement><textFieldExpression>$P{GreekText
} + "      " + $P{CyrillicText}</textFieldExpression></textField>
```

אני לא מבין עברית