



```
<image scaleImage="Clip" onErrorType="Icon" isLazy="true"><reportElement x="0"
y="0" width="150" height="40"/><imageExpression
class="java.lang.String">"http://jasperreports.sourceforge.net/jasperreports.png"</image
```



```
<image scaleImage="RetainShape"><reportElement x="0" y="50" width="150"
height="40"/><box
border="Thin"/><imageExpression>"dukesign.jpg"</imageExpression></image>
```



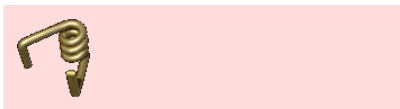
```
<image scaleImage="Clip"><reportElement x="0" y="100" width="150" height="40"
mode="Opaque" bgcolor="yellow"/><box border="Thin"
padding="5"/><imageExpression>"dukesign.jpg"</imageExpression></image>
```



```
<image scaleImage="FillFrame" isUsingCache="true"><reportElement x="0" y="150"
width="150" height="40"/><box
border="Thin"/><imageExpression>"dukesign.jpg"</imageExpression></image>
```



```
<image scaleImage="RetainShape" hAlign="Center"
isUsingCache="false"><reportElement x="0" y="200" width="150" height="40"/><box
border="Thin"/><imageExpression>"dukesign.jpg"</imageExpression></image>
```



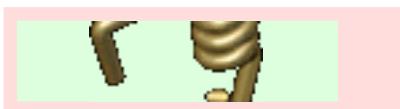
```
<image scaleImage="RetainShape"><reportElement x="5" y="255" width="70"
height="30"/><graphicElement/><imageExpression>"transparent.gif"</imageExpression
></image>
```



```
<image scaleImage="Clip"><reportElement x="5" y="305" width="70"
height="30"/><graphicElement/><imageExpression>"transparent.gif"</imageExpression
></image>
```



```
<image scaleImage="RetainShape"><reportElement x="5" y="355" width="30"
height="30" mode="Opaque"
bgcolor="#DDFFDD"/><graphicElement/><imageExpression>"transparent.gif"</imag
```



```
<image scaleImage="Clip" hAlign="Center" vAlign="Middle"><reportElement x="5"
y="405" width="120" height="30" mode="Opaque"
bgcolor="#DDFFDD"/><graphicElement/><imageExpression>"transparent.gif"</imag
```



```
<image scaleImage="Clip"><reportElement x="50" y="455" width="50" height="30"
forecolor="red"/><box
border="Thin"/><imageExpression>"transparent.gif"</imageExpression></image>
```



```
<image scaleImage="Clip"><reportElement x="50" y="505" width="50" height="30"
forecolor="red"/><box
border="Dotted"/><imageExpression>"transparent.gif"</imageExpression></image>
```



```
<image scaleImage="Clip"><reportElement x="50" y="555" width="50" height="30"
forecolor="red"/><box
border="1Point"/><imageExpression>"transparent.gif"</imageExpression></image>
```



```
<image scaleImage="Clip"><reportElement x="50" y="605" width="50" height="30"
forecolor="red"/><box
border="2Point"/><imageExpression>"transparent.gif"</imageExpression></image>
```



```
<image scaleImage="Clip" hAlign="Right" vAlign="Top"><reportElement x="50" y="655"
width="50" height="30" forecolor="red"/><box
border="4Point"/><imageExpression>"transparent.gif"</imageExpression></image>
```



```
<image><reportElement x="50" y="705" width="6" height="6" forecolor="green"/><box
border="4Point"/><imageExpression>"1redpixel.gif"</imageExpression></image>
```