

The quick brown fox jumps over the lazy dog.	<staticText><reportElement x="0" y="50" width="150" height="40"/><text>The quick brown fox jumps over the lazy dog.</text></staticText>
The quick brown fox jumps over the lazy dog.	<staticText><reportElement x="0" y="100" width="150" height="40"/><textElement></textElement><text>The quick brown fox jumps over the lazy dog.</text></staticText>
The quick brown fox jumps over the lazy dog.	<staticText><reportElement x="0" y="150" width="150" height="40"/><textElement></textElement><text>The quick brown fox jumps over the lazy dog.
<i>The quick brown fox jumps over the lazy dog.</i>	<staticText><reportElement x="0" y="200" width="150" height="40"/><textElement></textElement><text>The quick brown
The quick brown fox jumps over the lazy dog.	<staticText><reportElement x="0" y="250" width="150" height="40"/><textElement></textElement><text>The quick brown fox jumps over the
The quick brown fox jumps over the lazy dog.	<staticText><reportElement x="0" y="300" width="150" height="40"/><textElement></textElement><text>The quick brown fox jumps over
<u>The quick brown fox jumps over the lazy</u>	<staticText><reportElement x="0" y="350" width="150" height="40"/><textElement></textElement><text>The quick
The quick brown fox jumps over the lazy	<staticText><reportElement x="0" y="400" width="150" height="40"/><textElement></textElement><text>The quick
The quick brown fox jumps over the lazy	<staticText><reportElement x="0" y="450" width="150" height="40" forecolor="red"/><textElement></textElement><text>The quick brown fox jumps over the lazy dog.</text></staticText>
The quick brown fox jumps over the lazy dog.	<staticText><reportElement x="0" y="500" width="150" height="40" forecolor="green" backcolor="#FFFF00" mode="Opaque"/><textElement></textElement><text>The quick
The quick brown fox jumps over the lazy dog.	<staticText><reportElement x="0" y="550" width="150" height="90" forecolor="blue" backcolor="#FFDD99" mode="Opaque"/><textElement textAlignment="Center" verticalAlignment="Middle"></textElement><text>The quick brown fox jumps over the lazy dog.</text></staticText>
The quick brown fox jumps over the lazy dog.	<staticText><reportElement x="0" y="650" width="150" height="90" forecolor="red" backcolor="#99DDFF" mode="Opaque"/><textElement textAlignment="Right" verticalAlignment="Bottom"></textElement><text>The quick brown fox jumps over the lazy dog.</text></staticText>