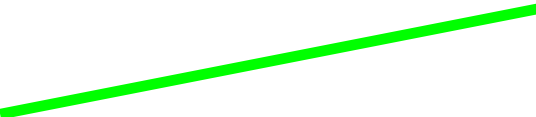




<line><reportElement x="0" y="50" width="200" height="0"/><graphicElement pen="1Point"/></line>



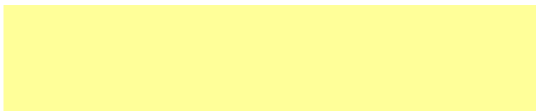
<line direction="TopDown"><reportElement x="0" y="100" width="200" height="40"/><graphicElement pen="2Point"/></line>



<line direction="BottomUp"><reportElement x="0" y="150" width="200" height="40" forecolor="green"/><graphicElement pen="4Point"/></line>



<line><reportElement x="100" y="200" width="0" height="40" forecolor="blue"/><graphicElement pen="Dotted"/></line>



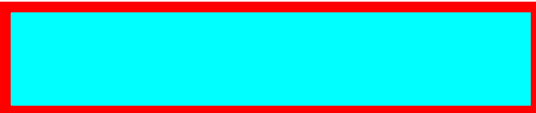
<rectangle><reportElement x="0" y="250" width="200" height="40" backcolor="#FFFF99"/><graphicElement pen="None"/></rectangle>



<rectangle><reportElement x="0" y="300" width="200" height="40"/><graphicElement pen="1Point"/></rectangle>



<rectangle radius="15"><reportElement x="0" y="350" width="200" height="40" backcolor="#FFFF00"/><graphicElement pen="2Point"/></rectangle>



<rectangle><reportElement x="0" y="400" width="200" height="40" forecolor="red" backcolor="#00FFFF"/><graphicElement pen="4Point"/></rectangle>



<rectangle><reportElement x="0" y="450" width="200" height="40" forecolor="#FF00FF" backcolor="#C0C0C0"/><graphicElement pen="Dotted"/></rectangle>



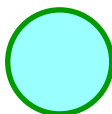
<rectangle><reportElement x="50" y="510" width="150" height="30" forecolor="blue" mode="Opaque"/><graphicElement pen="2Point"/></rectangle>



<rectangle><reportElement x="50" y="560" width="150" height="30" forecolor="blue" mode="Transparent"/><graphicElement pen="2Point"/></rectangle>



<ellipse><reportElement x="0" y="600" width="200" height="40" forecolor="red" backcolor="#FFFF99"/><graphicElement pen="2Point"/></ellipse>



<ellipse><reportElement x="80" y="650" width="40" height="40" forecolor="#009900" backcolor="#99FFFF"/><graphicElement pen="2Point"/></ellipse>