

100% shaded



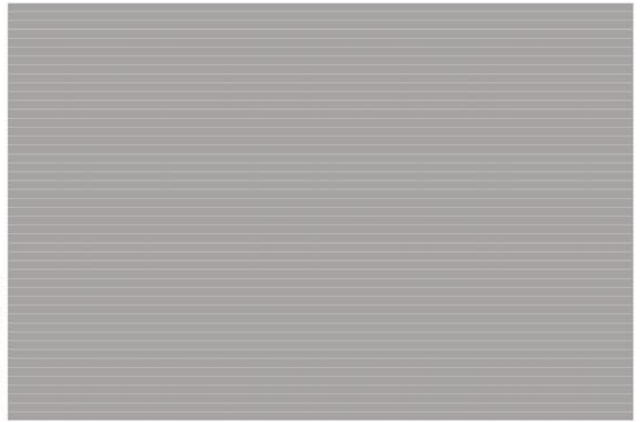
85% shaded



0% shaded



50% shaded



Red



Yellow



Green



Blue

