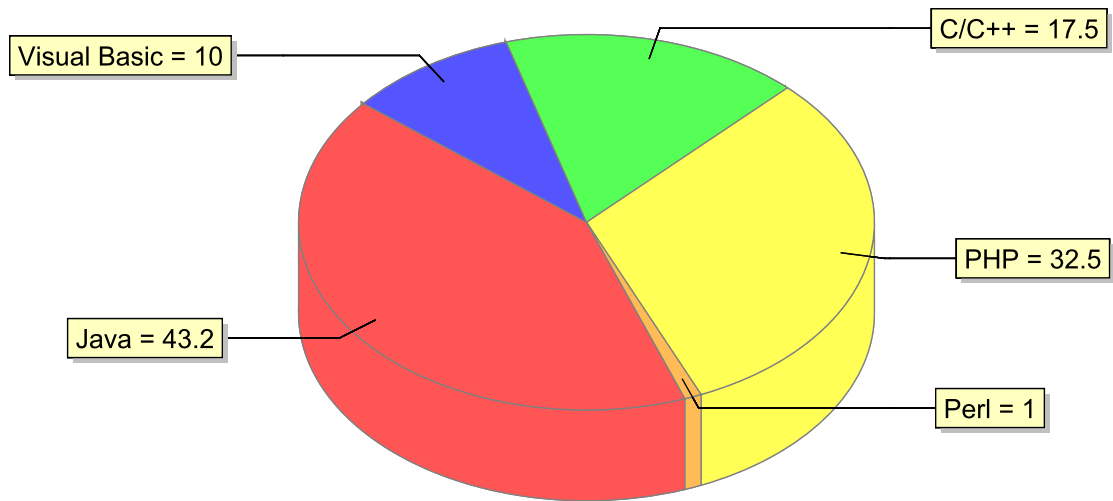


JFreeChart Sample

This sample uses JFreeChart Version 1.0.0-pre2
Written by David Gilbert (david.gilbert@object-refinery.com) and others.
(C)opyright 2000-2004, by Object Refinery Limited and Contributors.

Pie Chart 3D Demo 1



● Java = 43.2 ● Visual Basic = 10 ● C/C++ = 17.5 ● PHP = 32.5 ● Perl = 1