

---

# Barbecue Sample

This sample uses Barbecue Version 1.0.6b

<http://barbecue.sourceforge.net/>



1234567890