

DISCOVERING RICHARDIII

Turi King, University of Leicester

When the University of Leicester Archaeology Service undertook the Grey Friars project, it was thought that the chances of finding the remains of Richard III were slim to none.

Nevertheless, Turi King, with her background both in archaeology and genetics, was approached in the very early stages - should the skeletal remains of a 'good candidate' to be Richard III be found, would she be interested in overseeing the DNA analysis. On February 4th, 2013, Turi King was part of the team which presented the results of the analysis of Grey Friars Skeleton 1. In this talk, Dr Turi King will give an overview of the Grey Friars project.