

COMPARISON OF THREE DIFFERENT BLOOD EXTRACTION METHODS

K Tjhen
Bio-Synthesis, Inc.

This paper is to find out the advantages of three different blood extraction methods. These three methods are Chelex, FTA and whole blood extractions. The study is to extract the same blood sample with three different methods, and compare the data from electropherograms.