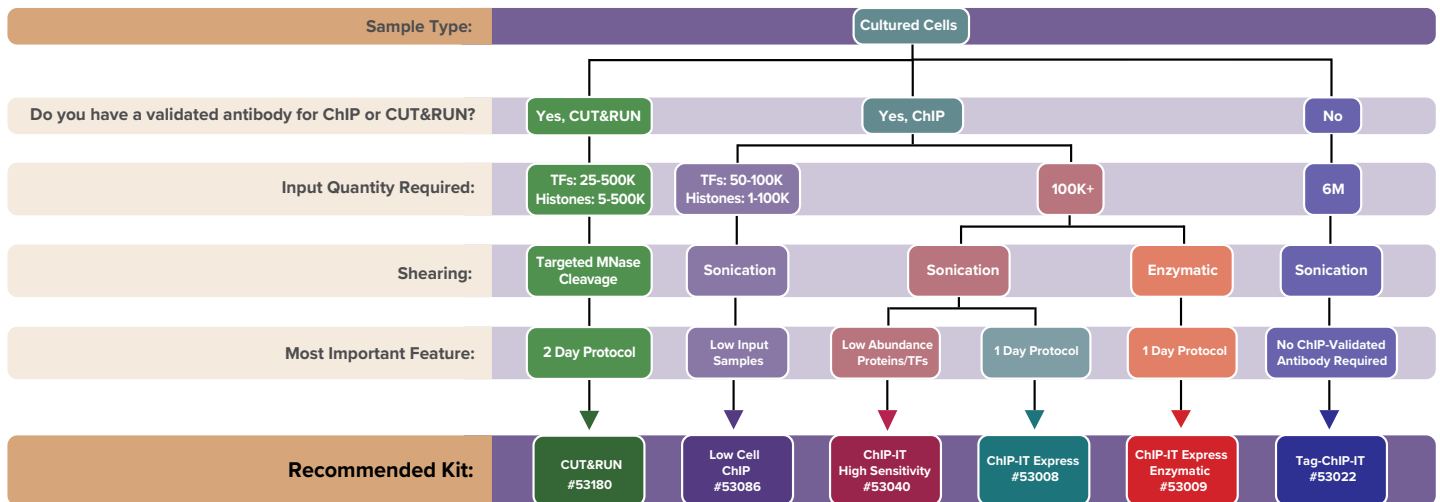


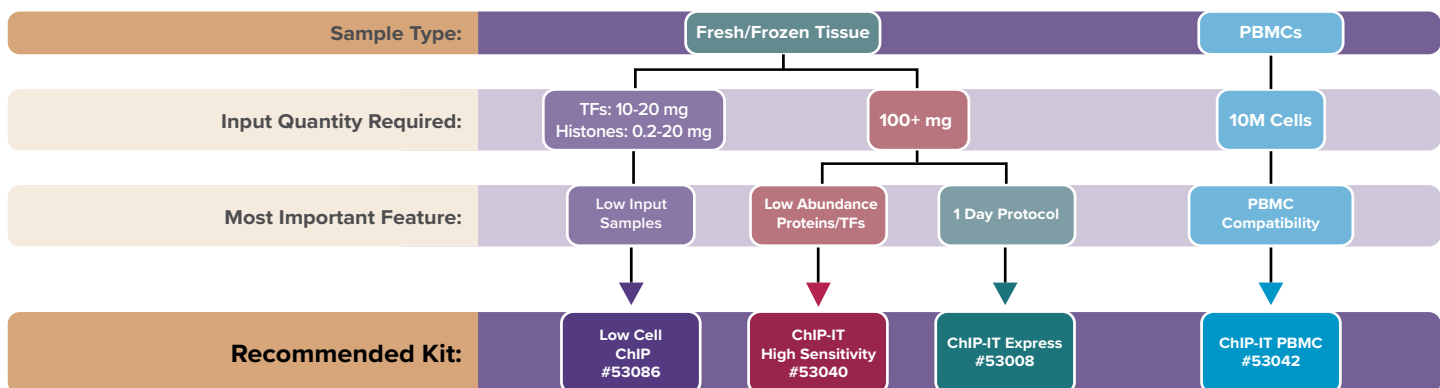
CHIP KIT SELECTION GUIDE

What is your starting material?

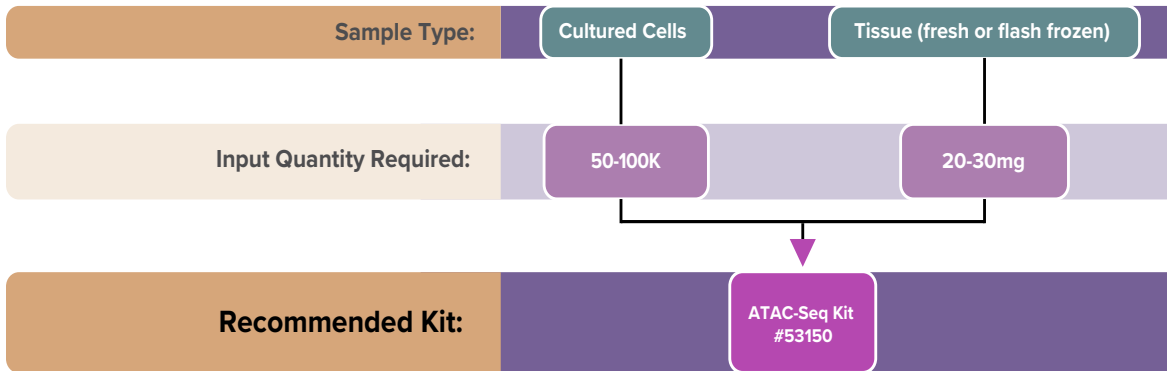
Cultured Cells



Fresh/frozen tissue or FFPE / PBMCs



Profile Open Chromatin Regions



Limited Sample Material

