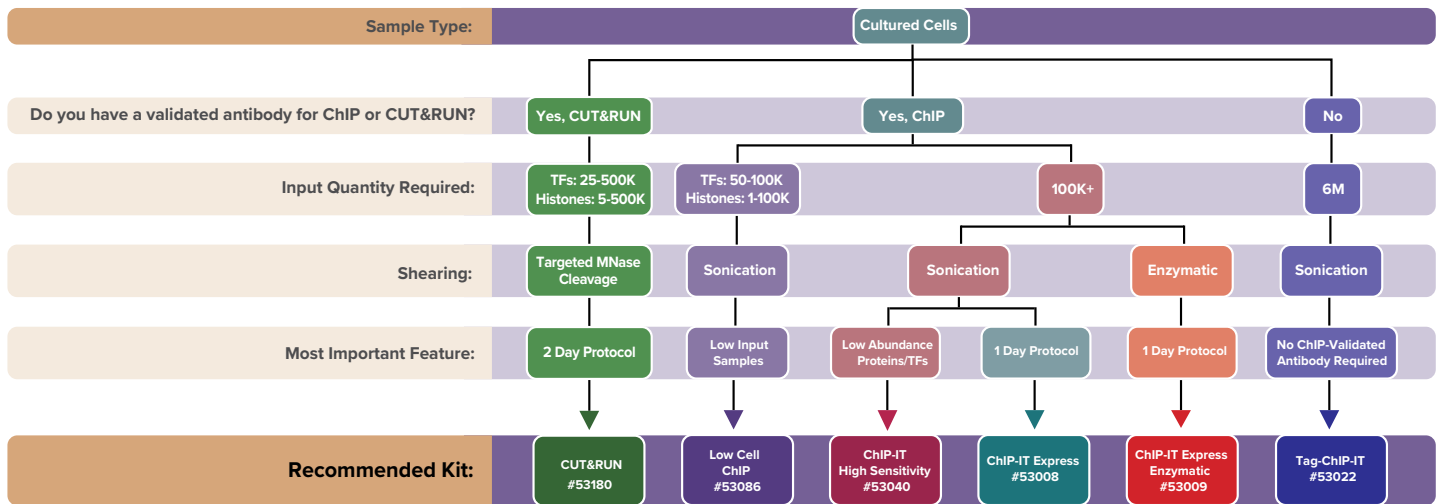


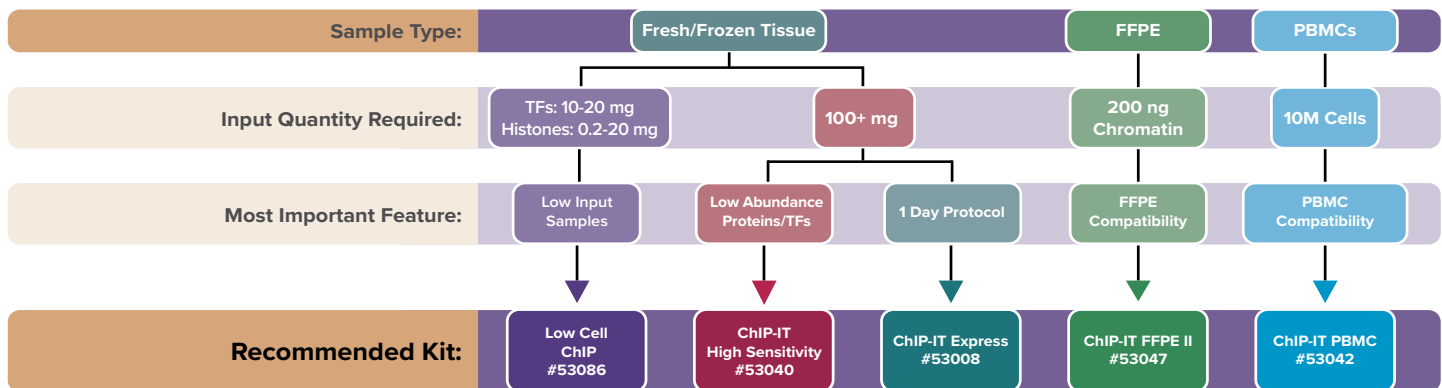
# CHIP KIT SELECTION GUIDE

## What is your starting material?

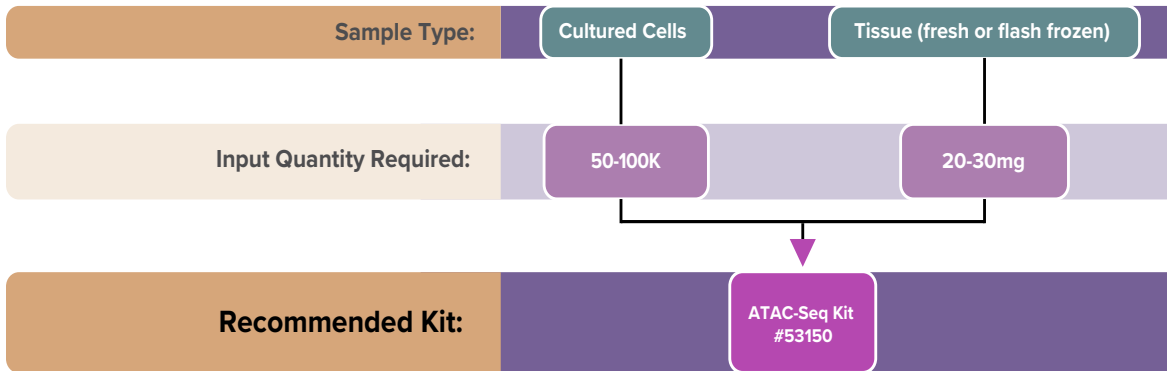
### Cultured Cells



### Fresh/frozen tissue or FFPE / PBMCs



## Profile Open Chromatin Regions



## Limited Sample Material

