

File : C:\ISO\05DEC2016\20803-0494659.D
Operator : SteveA
Acquired : 5 Dec 2016 19:45 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: 5-fluoro MDMB-PICA
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@240C,30C/min to 300C
Vial Number: 6

