

File :C:\ISO\12JAN2017\14764-0496720.D
Operator : SteveA
Acquired : 12 Jan 2017 19:21 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 5-fluoro ADB-PINACA
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 7

