

File :C:\ISO\16JAN2019 skg\9001123-0549363.D
Operator : SGoodwin
Acquired : 16 Jan 2019 13:11 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: bk-MDEA (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 4

