

File :C:\ISO\13MAY2016\9002651-0480761.D
Operator : SteveA
Acquired : 13 May 2016 15:42 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: (±)-threo-4-fluoromethylphenidate
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 11

