

File :C:\ISO\16MAY2017\14476-0504672R.D

Operator : JianmeiL

Acquired : 16 May 2017 13:41 using AcqMethod CAY_DRUG_CATH.M

Instrument : ISO_GCMS1

Sample Name: Sibutramine (HCl)

Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C

Vial Number: 5

