

File :C:\ISO\Data\30JUN2015\17735-0469229R.D
Operator : LauraN
Acquired : 1 Jul 2015 15:19 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 5-fluoro AEB
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 8

