

File : C:\ISO\01DEC2016\12063-0480732.D  
Operator : SteveA  
Acquired : 1 Dec 2016 15:11 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : ISO\_GCMS1  
Sample Name: etizolam  
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C  
Vial Number: 3

