

File : C:\ISO\30MAR2017\9001355-0459577r.D 0459577 -SKB17AUG2018  
Operator : SamanthaB  
Acquired : 31 Mar 2017 15:19 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : ISO\_GCMS1  
Sample Name: 3-methoxy PCE (hydrochloride)  
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C  
Vial Number: 7

