

File :C:\ISO\20SEP2018 skb\15919-0540757 R.D  
Operator : SBouhall  
Acquired : 20 Sep 2018 13:12 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : ISO\_GCMS1  
Sample Name: cis-Tramadol (hydrochloride)  
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C  
Vial Number: 6

