

File :Y:\Forensics\9001008-0439311.D  
Operator : PDK  
Acquired : 10 Aug 2012 22:15 using AcqMethod CAY\_DRUG.M  
Instrument : Instrument #1  
Sample Name: JWH 018 2'-naphthyl-N-(2,2-dimethylpropyl) is  
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C  
Vial Number: 14

